

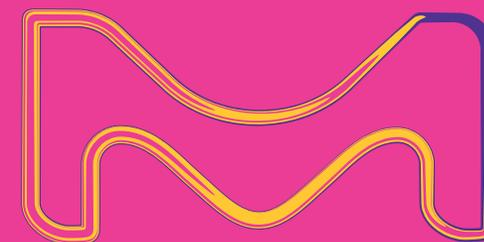
**SAFC®**

Pharma & Biopharma  
Raw Material Solutions

# Strategies and Enabling Technologies for Enhancing API Solubility

Dr. Markus Lubda, Strategic Marketing Manager  
Dr. Thomas Briel, Strategic Marketing Manager

Webinar – April 2023



The Life Science business  
of Merck KGaA, Darmstadt, Germany operates as  
MilliporeSigma in the U.S. and Canada



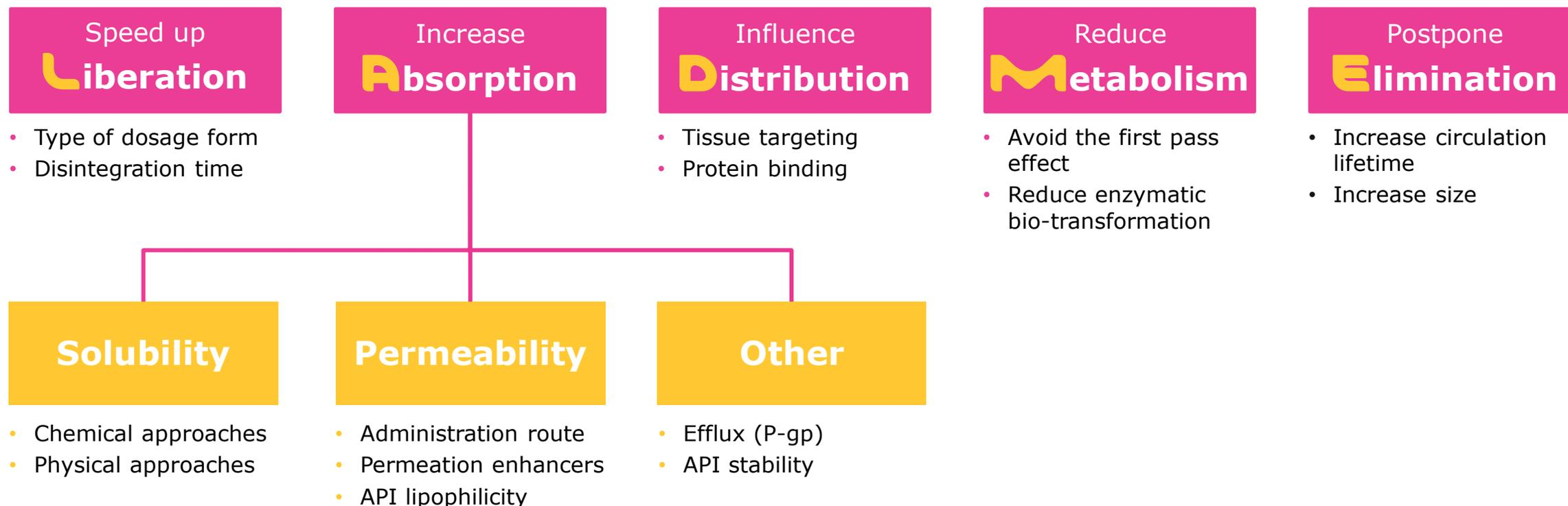
# Agenda

- 01 Introduction to Developability Classification System (DCS)
- 02 Strategies to enhance solubility of DCS Class IIa molecules
- 03 Strategies to enhance solubility of DCS Class IIb molecules
- 04 Summary



## Drug exposure

# The exposure of the drug in the bloodstream, is essential for a physiological effect

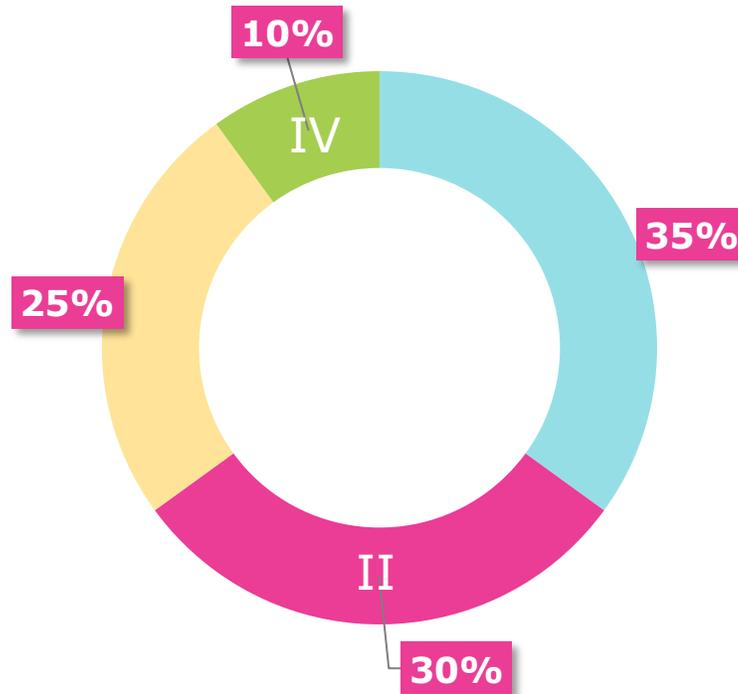
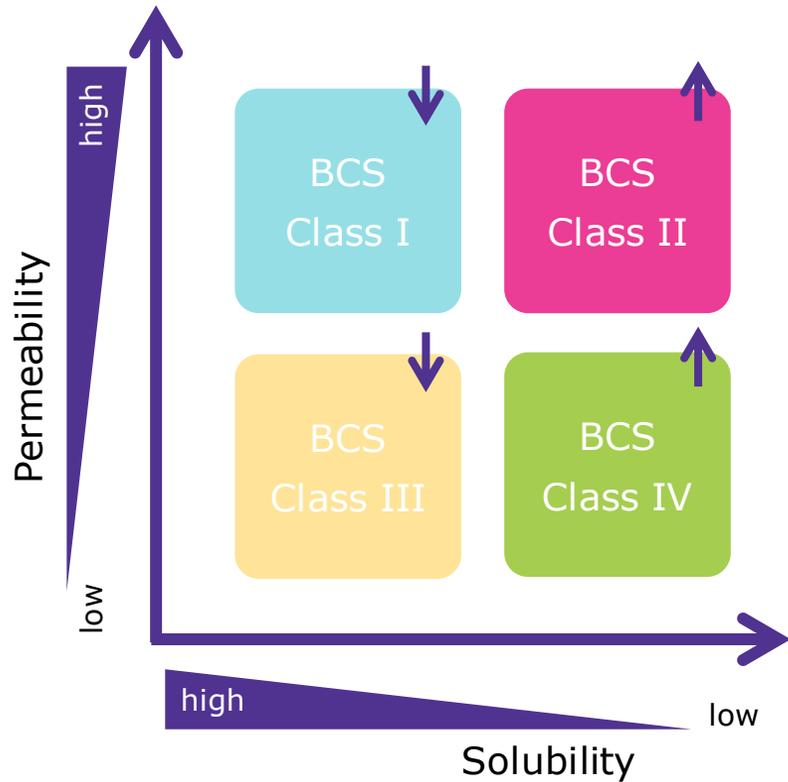


Absorption of the API from the intestinal tract to systemic circulation is governed primarily by **solubility and permeability**

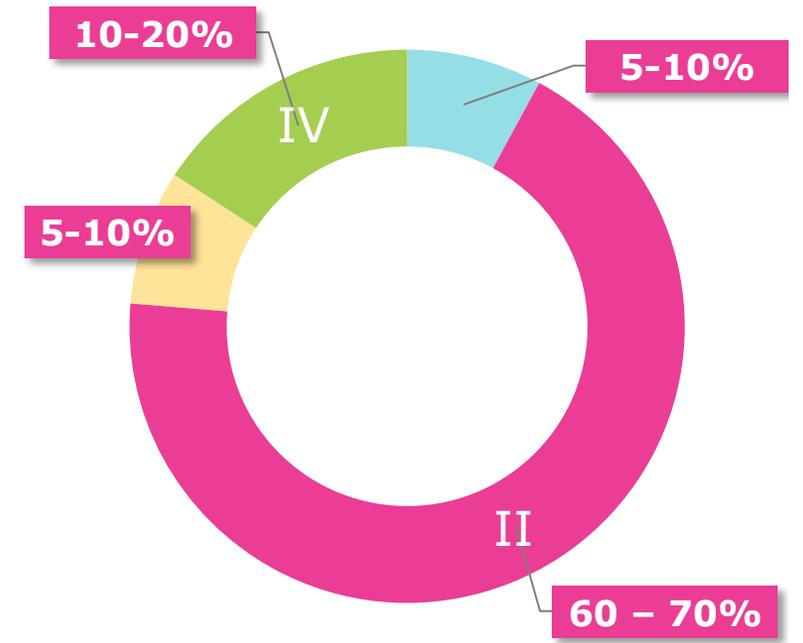


# Solubility of APIs in development

## Poor soluble molecules are becoming more prevalent



Distribution of **marketed** drug substances



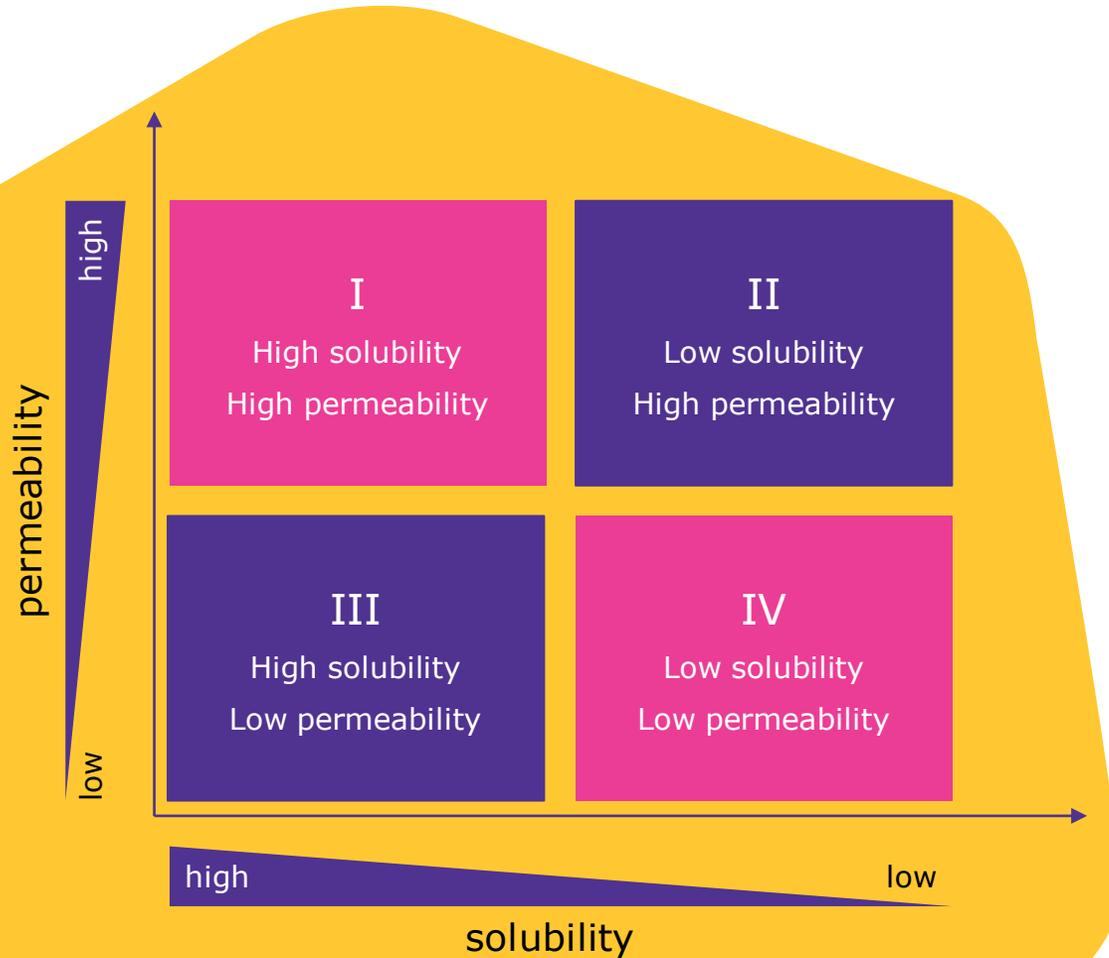
Distribution of **pipeline** drug substances



**Poor solubility can lead to low and variable absorption**



## The BCS classifies molecules based on solubility and permeability



### Originated in the 1990s

Correlates *in-vitro* to *in-vivo* performance based on aqueous solubility and permeability, and classifies APIs based on these factors

### Adopted by regulatory Authorities

Used by FDA as basis for so called „Biowaiver“ applications for immediate-release formulations in order to reduce unnecessary additional in-vivo studies for bioavailability and bioequivalence.

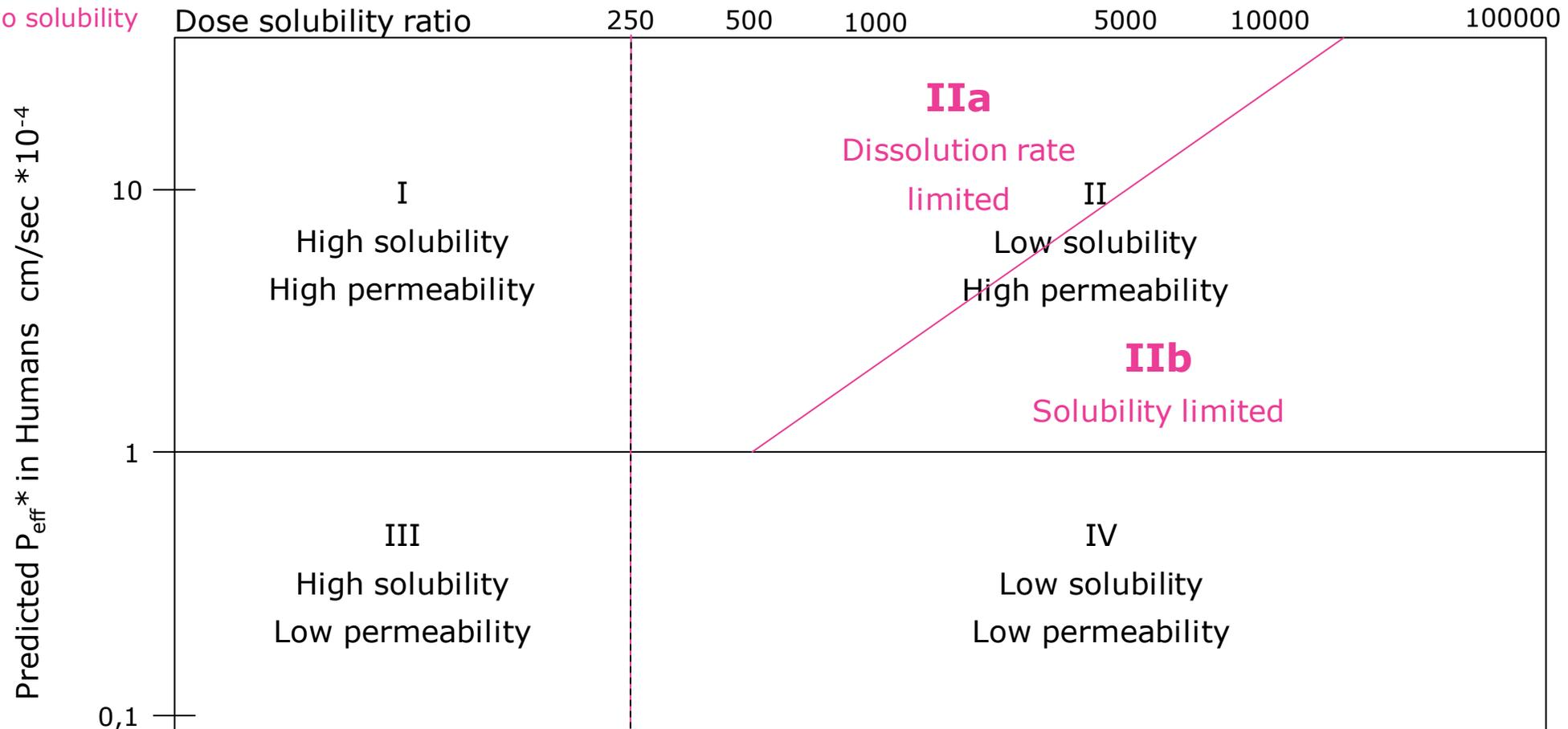


# The Developability Classification System (DCS)

## The DCS is based on the BCS and can be used to optimize formulations of poorly soluble molecules

Using FaSSIF for an estimate

of in vivo solubility



\* Effective intestinal membrane permeability



# General approach

## Applying DCS in formulation development

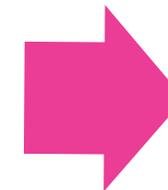
### Solubility Determination

- Thermodynamic solubility
- Describes the maximum saturation concentration
- Measured in FaSSIF
- Described using the Dose/Solubility ratio, the minimum amount of “solvent” required to dissolve the maximum dose



### Permeability Determination

- Description of flux through cell membrane
- Measured using Caco-2 cells
- Can also be predicted using software



### DCS Class Assignment

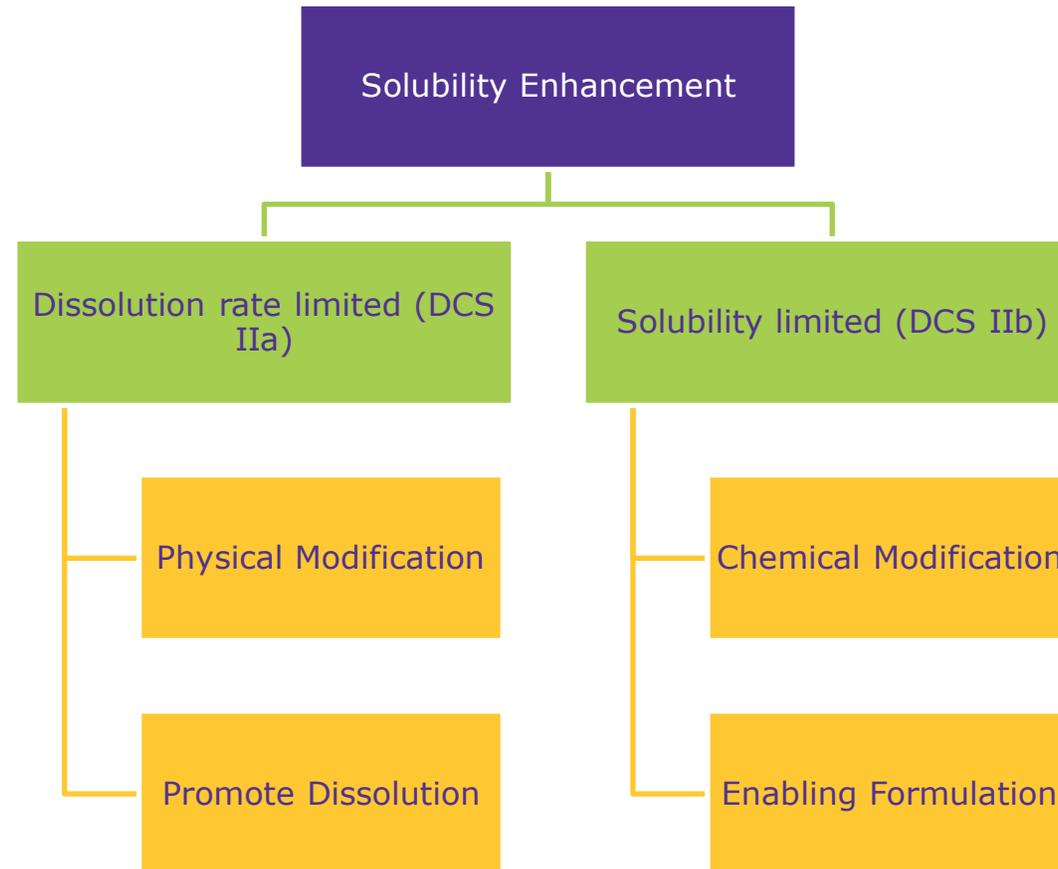
- Based on the combination of solubility and permeability
- Determines its DCS class
- For DCS II molecule, divided by the solubility limited absorbable dose (SLAD) line



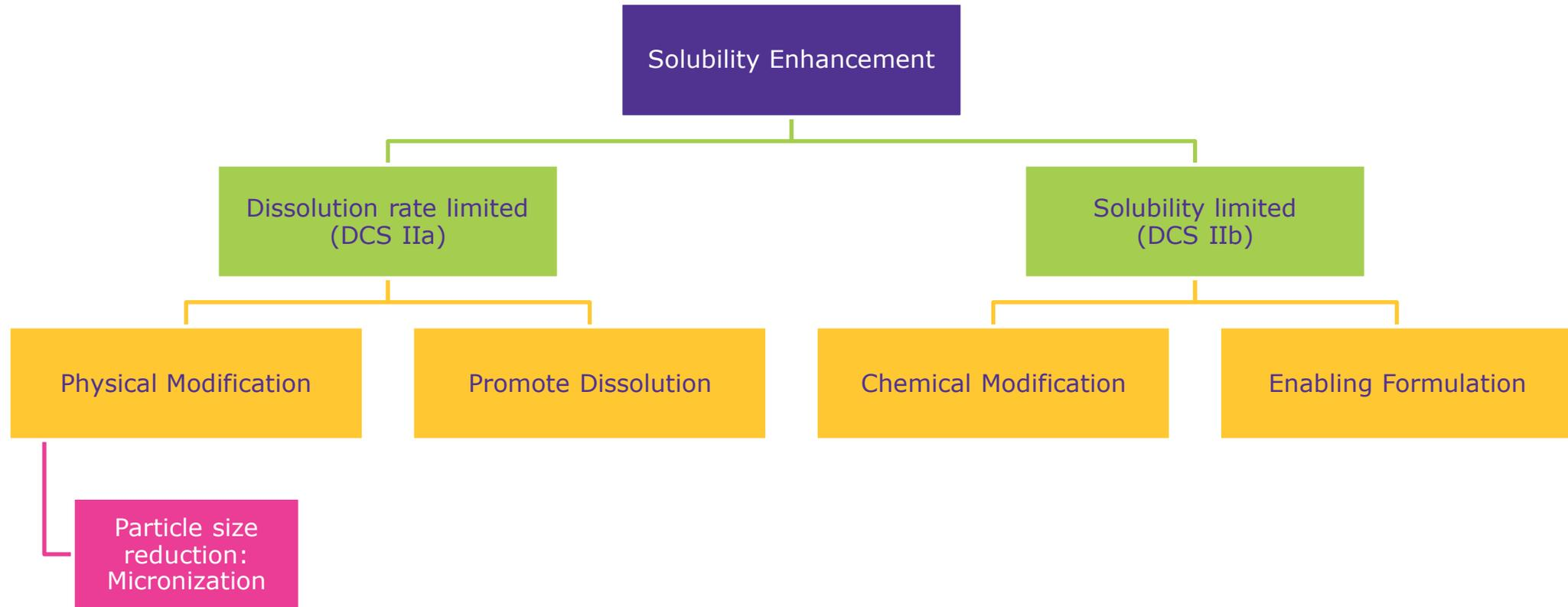
**The DCS class of a molecule can be determined using readily accessible laboratory methods and equipment**



# Strategies and Enabling Technologies for Enhancing API Solubility



# Strategies and Enabling Technologies for Enhancing API Solubility



- According to Noyes-Whitney equation, dissolution rate of a drug particle is related to its surface area  
→ **smaller particles means higher dissolution**<sup>[1]</sup>
- Micronization aims at reducing particle size to  $\mu\text{m}$  range, but no generally defined borders
- Frequently used methods are **mechanical milling** (coarse particles) and **jet milling** (fine particles)<sup>[2]</sup>

## Mechanical Milling

- Typically particle sizes in the range 50 – 75  $\mu\text{m}$
- Established, easy to handle, rather cheap
- Moving / rotating parts with mechanical wear
- Abrasion of machinery possible -> contamination
- Limited control of particle size distribution

## Jet Milling

- Particle sizes down to approx. 1  $\mu\text{m}$  possible
- Employs pressurized gas at >sonic speed
- No moving parts, low / no mechanical wear
- Contamination risk strongly reduced
- Higher control of particle size distribution

[1] The rate of solution of solid substances in their own solutions, Noyes & Whitney, 1897

[2] Overview of nanoparticulate strategies for solubility enhancement of poorly soluble drugs, Khan *et al.*, 2022



## ***In-Situ* Micronization**

- Principle: Drug is precipitated with a non-solvent, followed by drying
- Poorly water-soluble drugs: solubilized with an organic solvent; aqueous solvent (including stabilizing agent) is used as precipitator
- Common stabilizers: PEG, PVA, PVP, Cellulose (API-dependent)
- Particle sizes down to approx. 1  $\mu\text{m}$  possible

## **Benefits:**

- No specialized mechanical equipment needed
- Tight particle size distribution, drug amount in preparation relatively high<sup>[2]</sup>
- Polymer-hydrophilized surface increases dissolution and prevents aggregation

## **Milling**

- Well-established techniques
- Thermal and physical stress
- Mechanical equipment needed
- Size control depends on method

## ***In-Situ* Micronization**

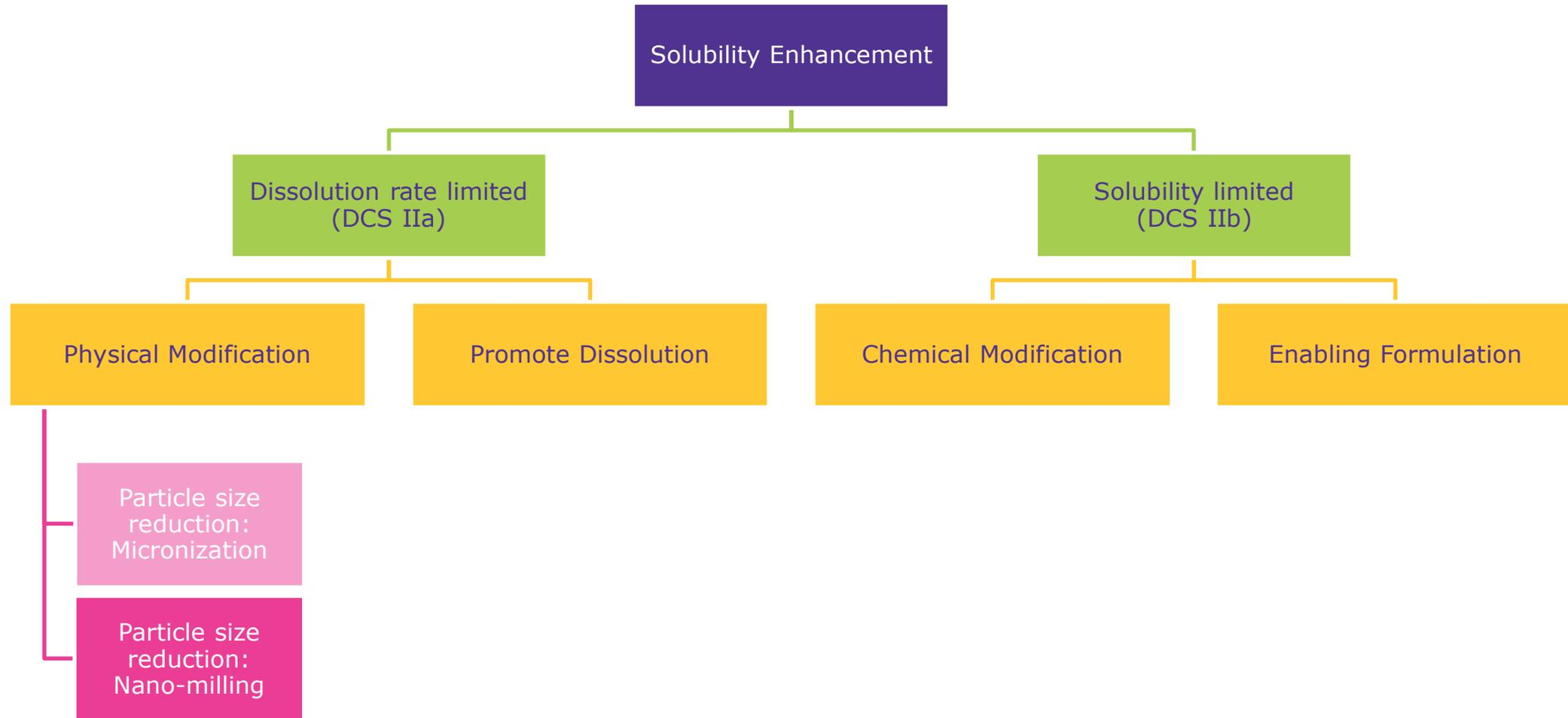
- No stress, tight size control
- Process rather demanding
- Stabilizers needed

[1] An overview on in situ micronization technique – An emerging novel concept in advanced drug delivery, Vandana *et al.*, 2013

[2] Dissolution rate enhancement by in situ micronization of poorly water-soluble drugs, Rasenack & Müller, 2002

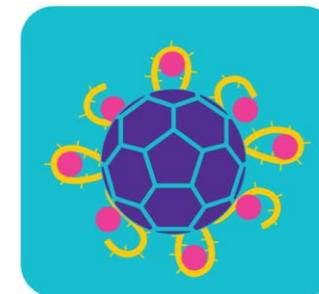
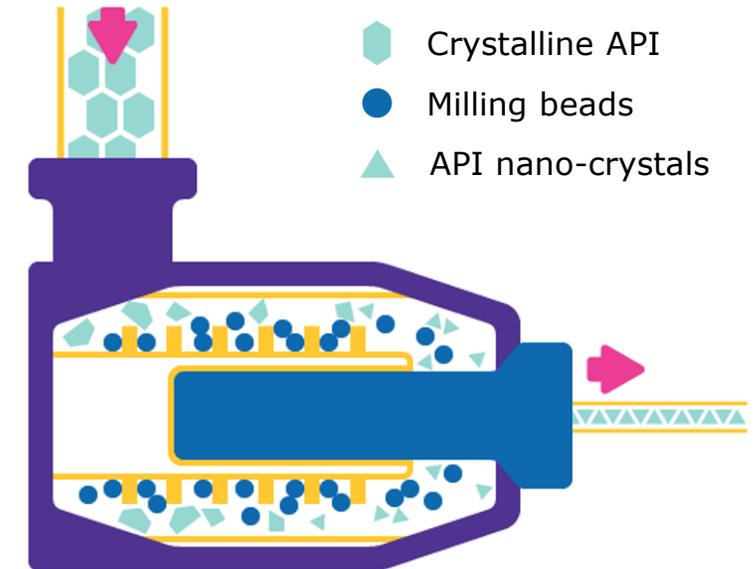


# Strategies and Enabling Technologies for Enhancing API Solubility



# Nano-Milling

- API **nano-milling is an established procedure** to enhance dissolution by strongly increasing the drug surface area<sup>[1]</sup>
- Target is usually a reduction of **particle size below 1000 nm**, typically 100 nm range
- **Can be converted into diverse dosage forms** like oral liquids, capsules, tablets, injectables, aerosols, etc.<sup>[1]</sup>
- Milling is **performed using a medium with beads** made of ceramics or highly crosslinked polystyrene<sup>[2]</sup>
- **Outcome is a nanosuspension** comprised of water, stabilizer(s), and API



stabilized drug  
nano-crystal

[1]: Nanomilling of Drugs for Bioavailability Enhancement: A Holistic Formulation-Process Perspective, Li *et al.*, 2016

[2]: Nanomilling: A Key Option for Formulating Water-Insoluble APIs, Lee & DiFranco, 2022



## Examples Marketed Nano-Milled Drugs<sup>[1]</sup>:

Product	API	Company
Rapamune <sup>®</sup>	Rapamycin, Sirolimus	Wyeth
Emend <sup>®</sup>	Aprepitant	Merck & Co.
Tricor <sup>®</sup>	Fenofibrate	Abbott Laboratories
Ritalin LA <sup>®</sup>	Methylphenidate HCl	Novartis
Zanaflex <sup>®</sup>	Tizanidine HCl	Acorda
Invega Sustenna <sup>®</sup> / Trinza <sup>®</sup>	Paliperidone Palmitate	Janssen Pharma
Yonza <sup>®</sup>	Abiraterone Acetate	Sun Pharma

## Due to high re-aggregation tendency stabilizers are required:

Poloxamer	Polyethylene glycol
Tween	Polyvinyl alcohol
Sodium dodecyl sulfate	

### Benefits of Nano-Milling

- Well established
- No chemical modification
- No limitation for dosage form

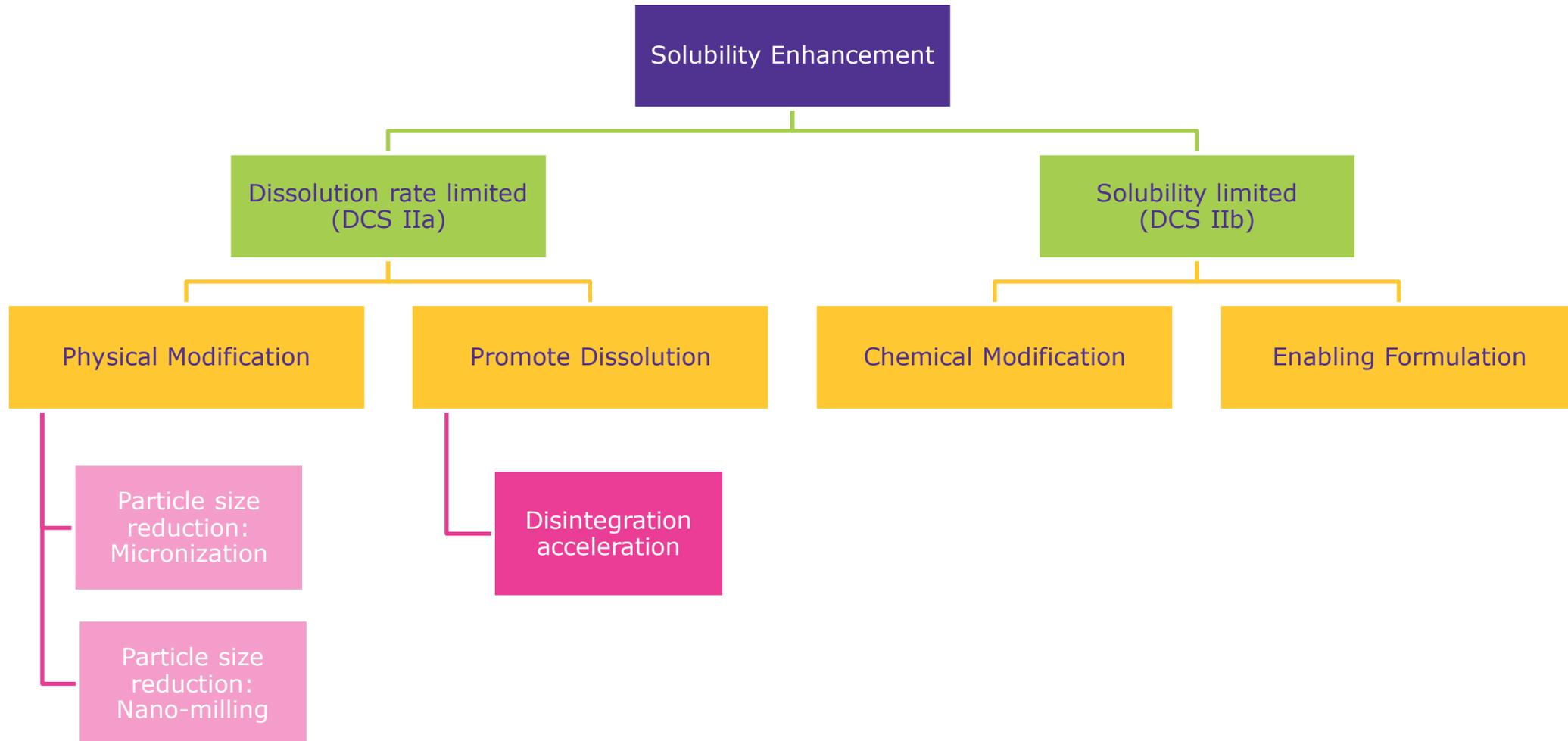
### Challenges

- Mechanical stress to API
- Re-aggregation can occur
- Stabilizers needed

[1]: Nanomilling: A Key Option for Formulating Water-Insoluble APIs, Lee & DiFranco, 2022



# Strategies and Enabling Technologies for Enhancing API Solubility



Accelerate disintegration

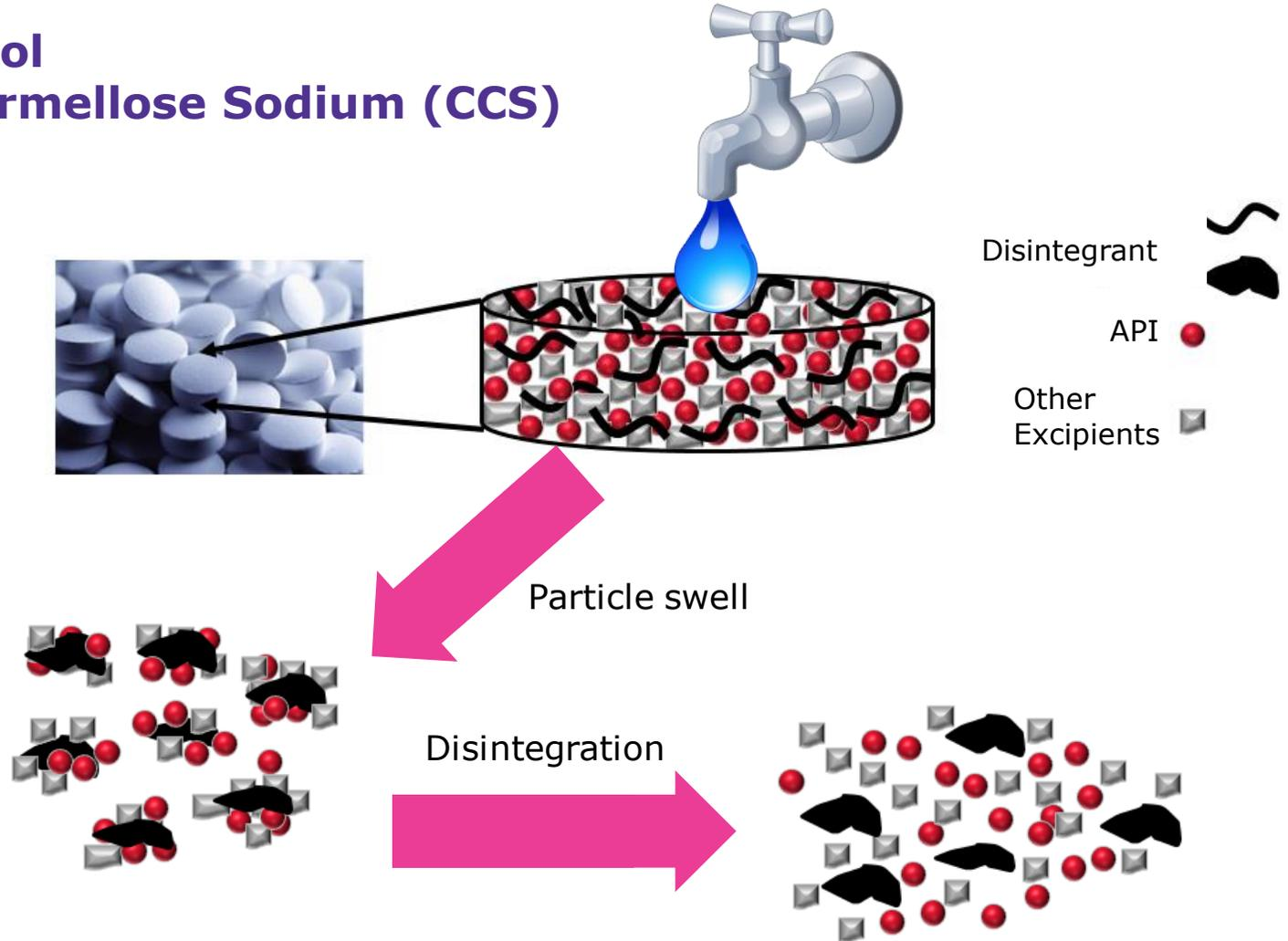
**Disintegrants are important for the dissolution rate**

➤ **Mannitol**  
**Croscarmellose Sodium (CCS)**

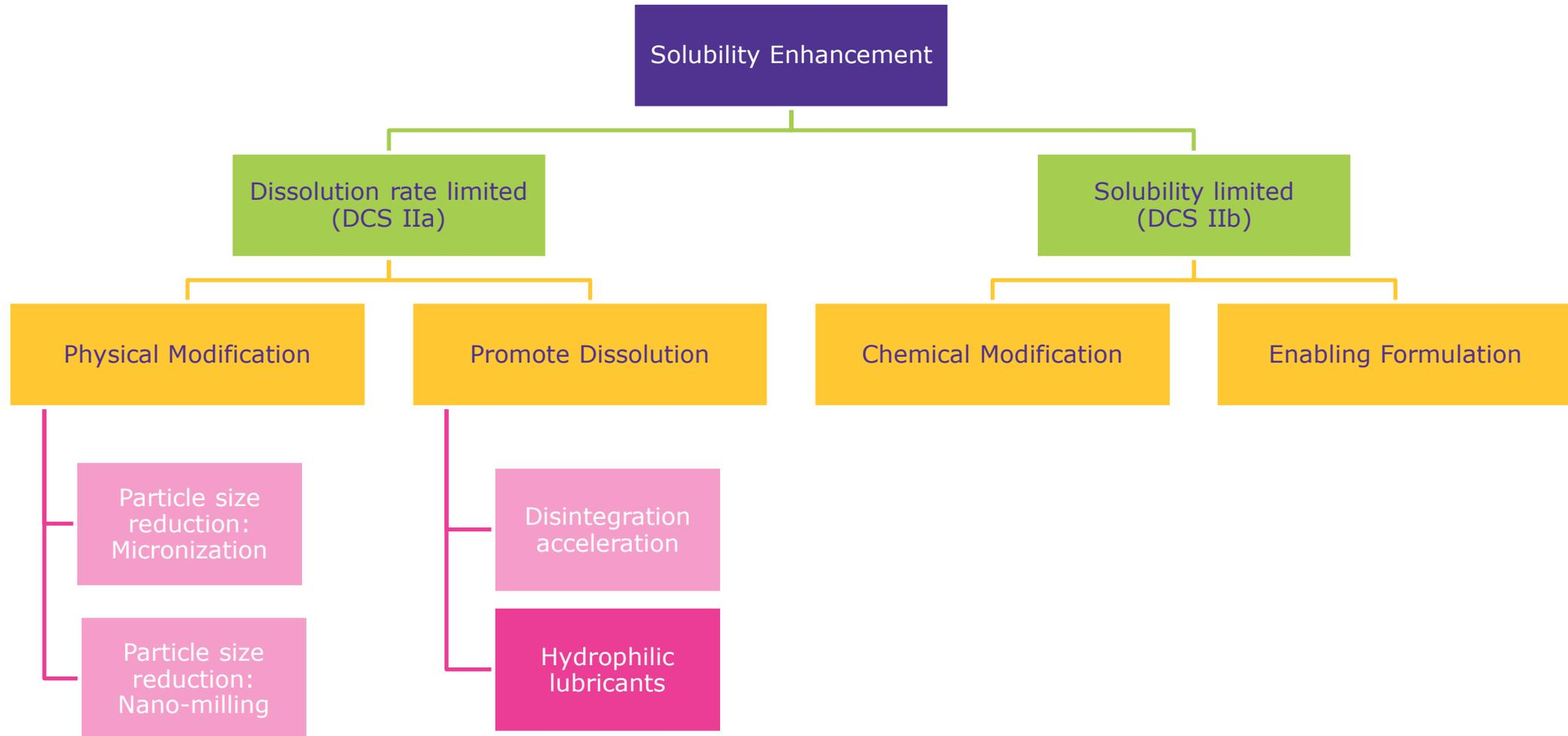
**Disintegrant features**

- Good hydration capacity
- Surface structure
- Compressibility
- Compatible with enabling technology

➔ **Fast disintegration**



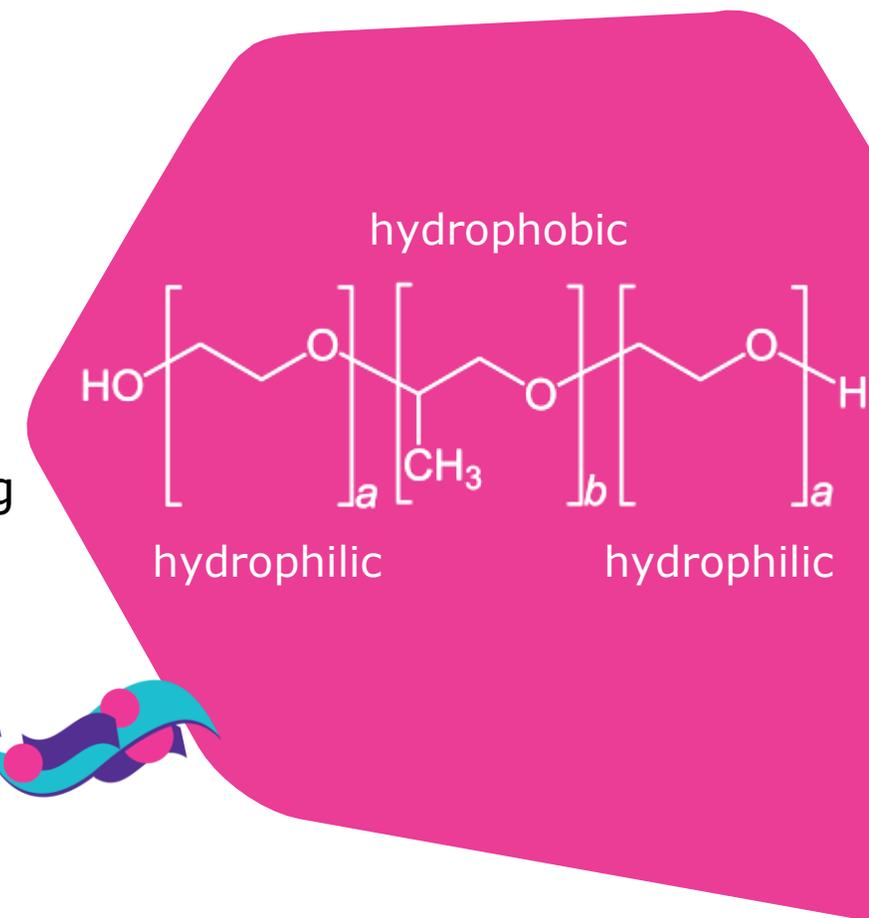
# Strategies and Enabling Technologies for Enhancing API Solubility



## Hydrophilic Lubricant – Dissolution enhancement

### A multi-functional excipient

- **Poloxamer 188**, a block copolymer of poly(ethylene oxide) and poly(propylene oxide)
- Functional excipient for oral solid dosage forms acts as a **dissolution rate enhancer** via a surfactant like effect
- Additional functionality as a **hydrophilic lubricant**
- Compatible for: DC, continuous manufacturing and 3D printing



Parteck® PLX 188

Dissolution enhancer 1 – 15 %

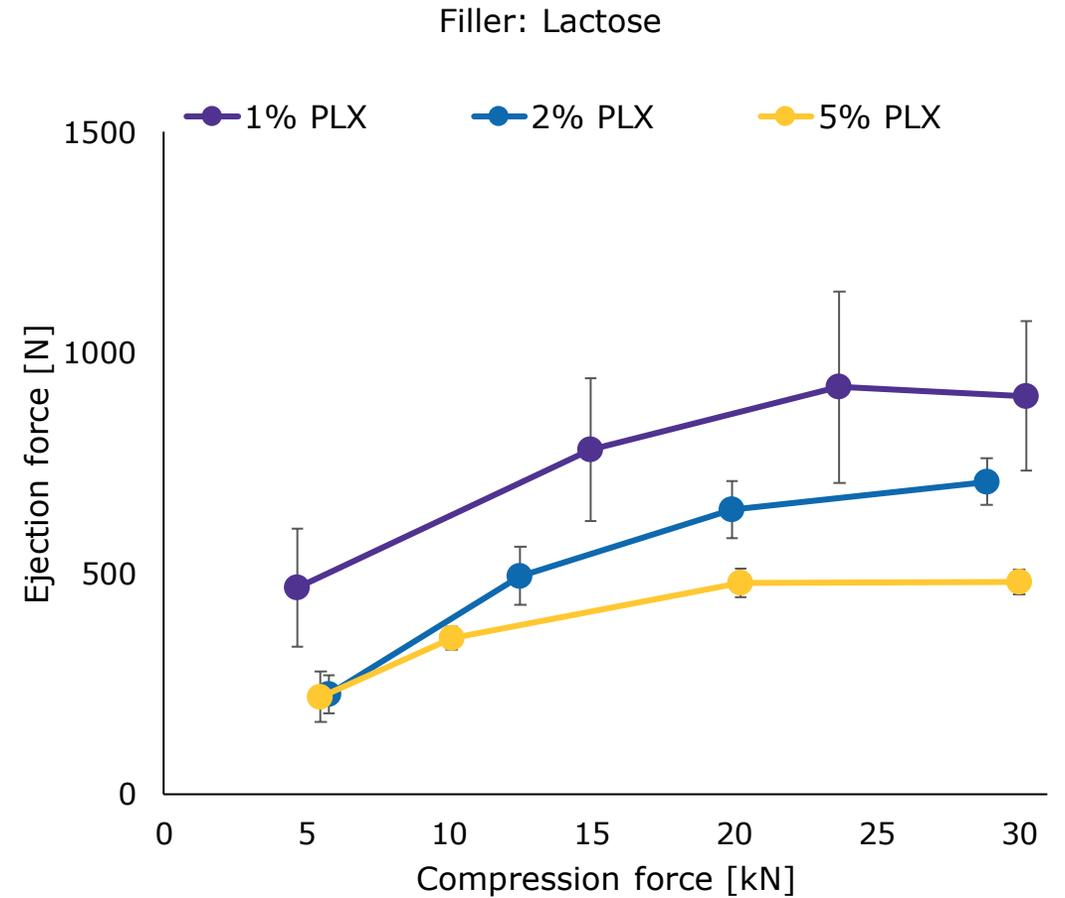
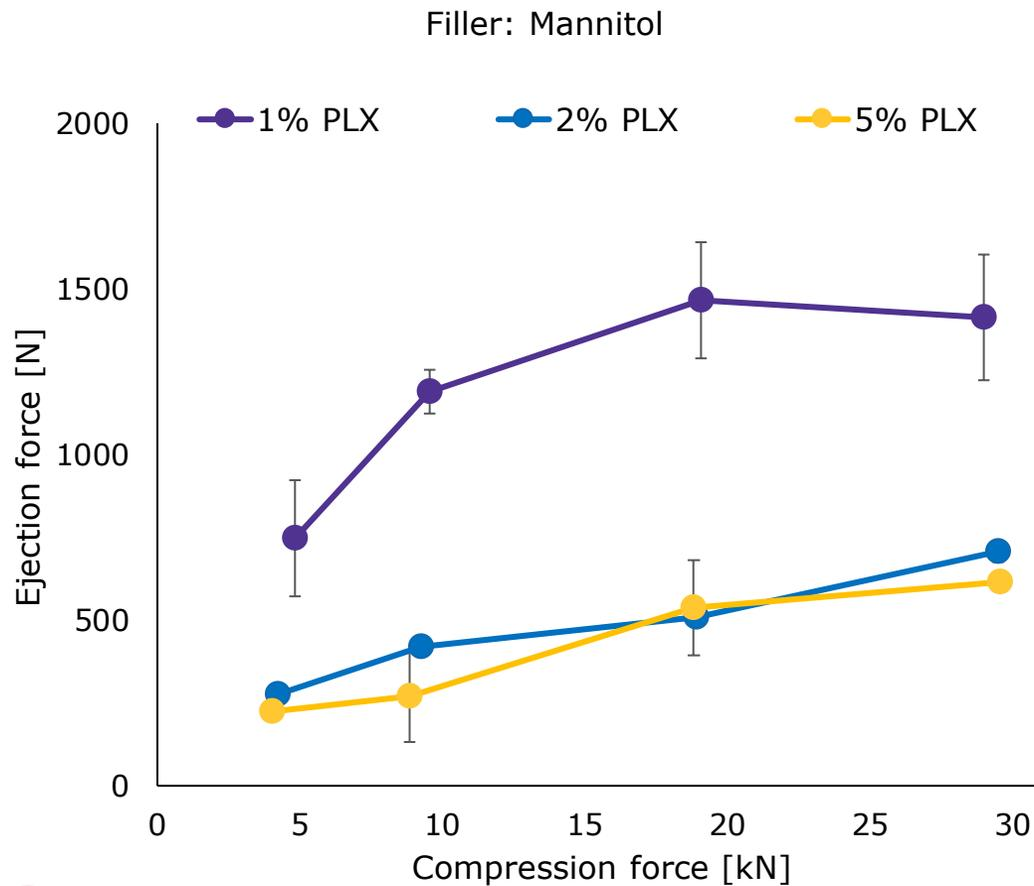
Lubricant 2 – 5 %

➤ **One excipient, multiple functionalities:**  
**Improving process efficiency and unlock future manufacturing capabilities**



# Poloxamer 188 – Hydrophilic lubricant

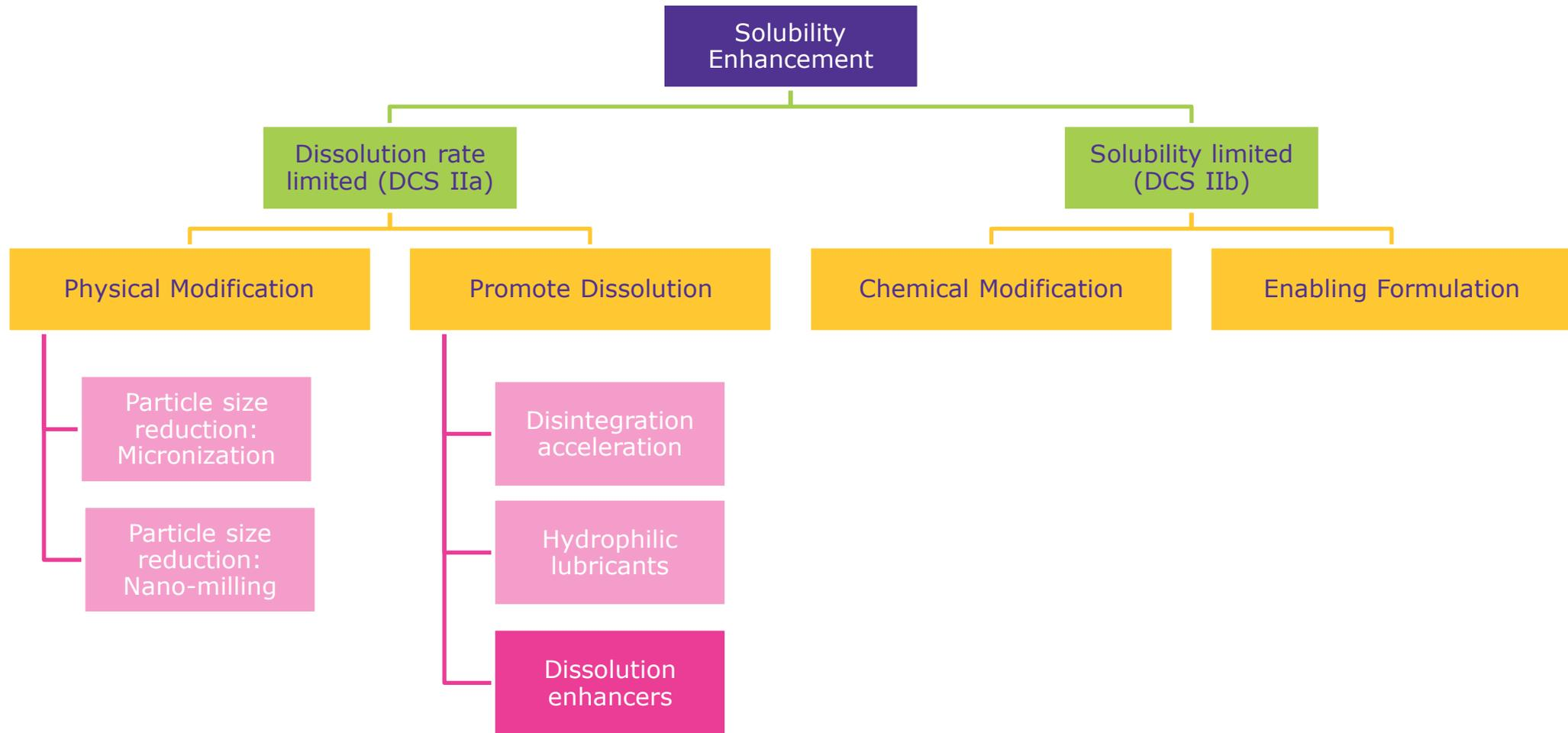
## Effective lubrication is achieved even when combined with brittle, deforming fillers



A minimum of 2% Poloxamer 188 is recommended for lubrication effect  
→ Perfect fit for direct compression

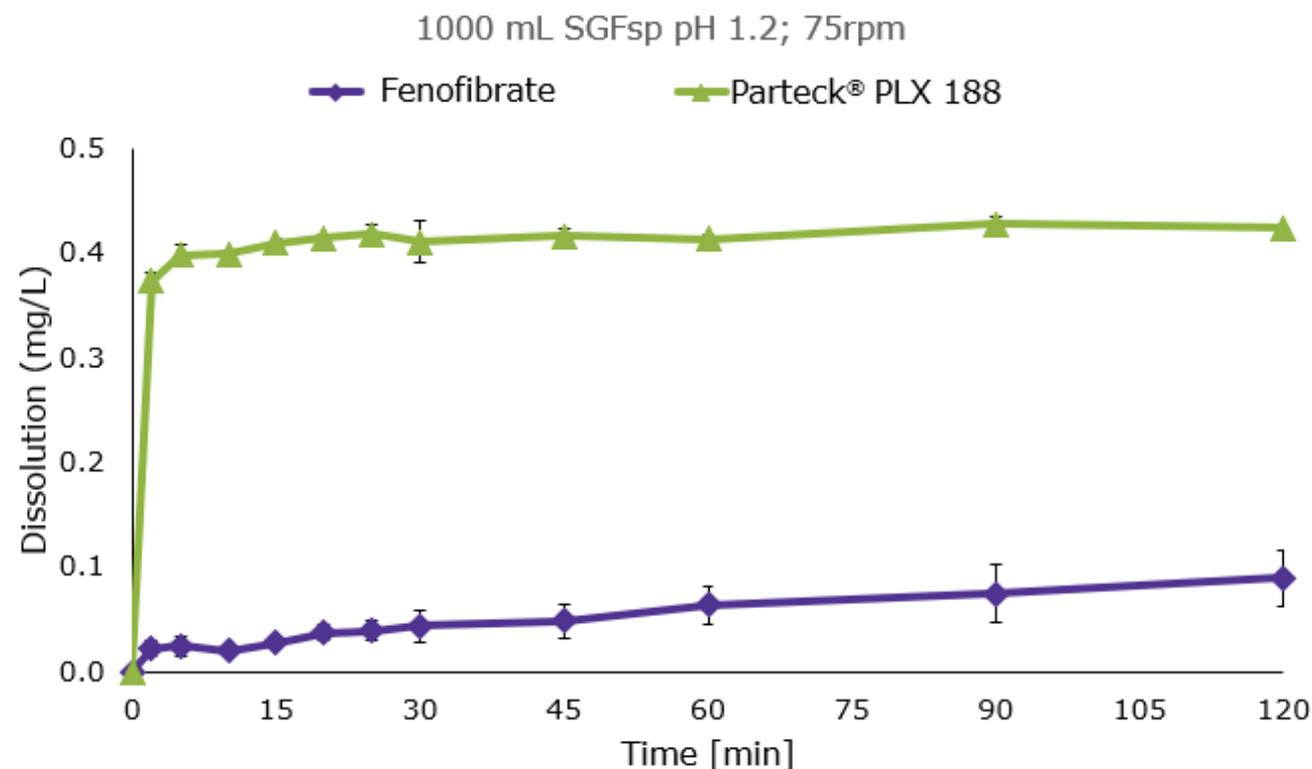


# Strategies and Enabling Technologies for Enhancing API Solubility



## Poloxamer 188 – Dissolution enhancement

### Dissolution is enhanced via increased wetting and micelle formation



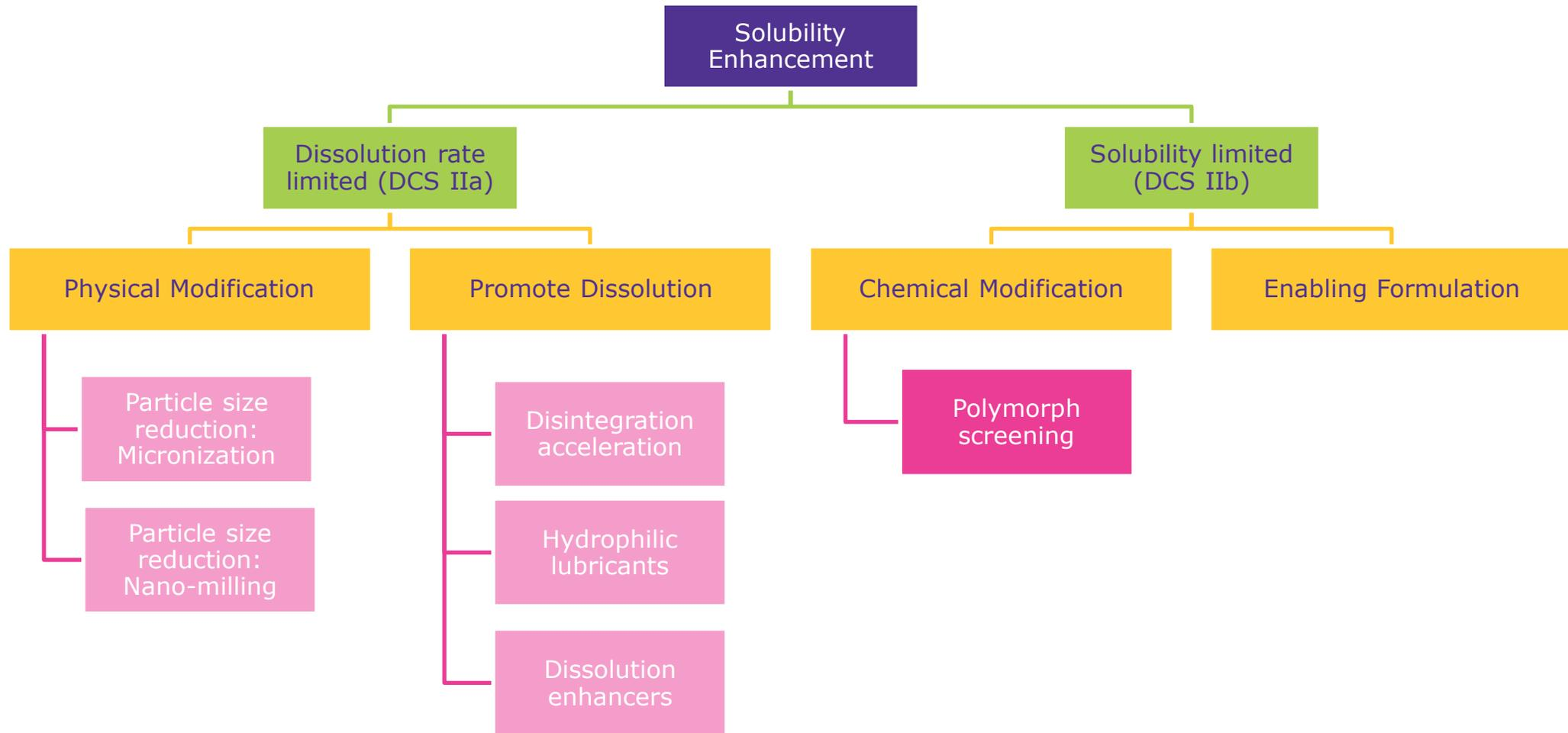
Component	Content [%]
<b>Parateck® PLX 188</b>	<b>5</b>
<b>Fenofibrate</b>	<b>20</b>
Parateck® M 200	69
Crosscarmellose sodium (CCS)	5
Magnesium stearate	1



**Adding Poloxamer 188 to the formulation is a fast and easy solution to enhance the dissolution rate**

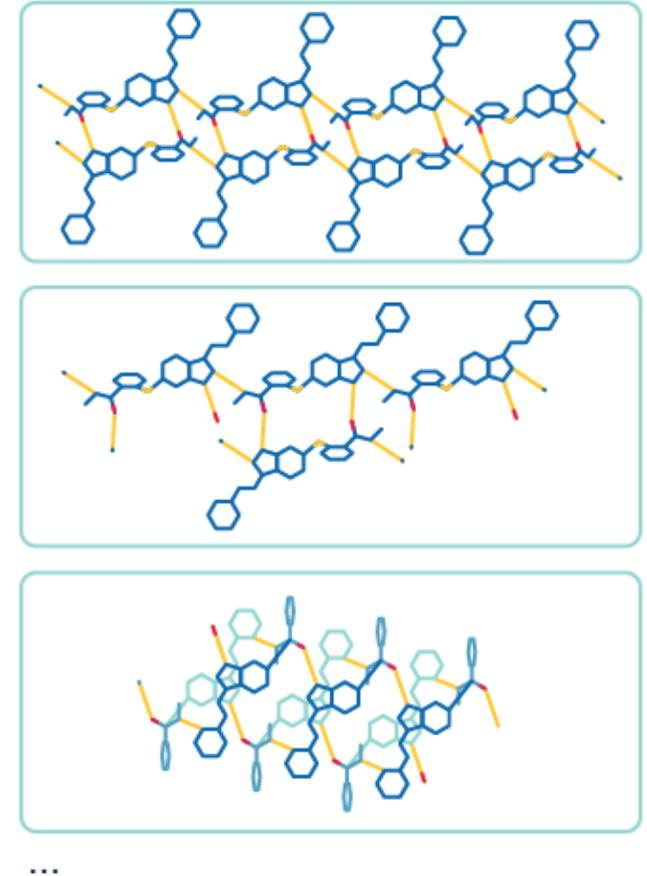


# Strategies and Enabling Technologies for Enhancing API Solubility



# Polymorph Screening

- **Polymorphism** means that drugs can have several crystalline forms – which is true for **>50 % of APIs**<sup>[1]</sup>
- Polymorph screening is **mandatory due to safety reasons**, e.g. regarding thermodynamic stability and toxicity (-> ICH Q6A)
- Polymorphs can be considerably different regarding<sup>[2]</sup>
  - Solubility
  - Biological activity
  - Melting Point
  - Dissolution rate
  - Pharmaco-dynamics
- Polymorphic forms are considered **different chemical entities**



[1] Development of a Targeted Polymorph Screening Approach for a Complex Polymorphic and Highly Solvating API, Campeta *et al.*, 2010

[2] Polymorphism: The Phenomenon Affecting the Performance of Drugs, Raza *et al.*, 2014

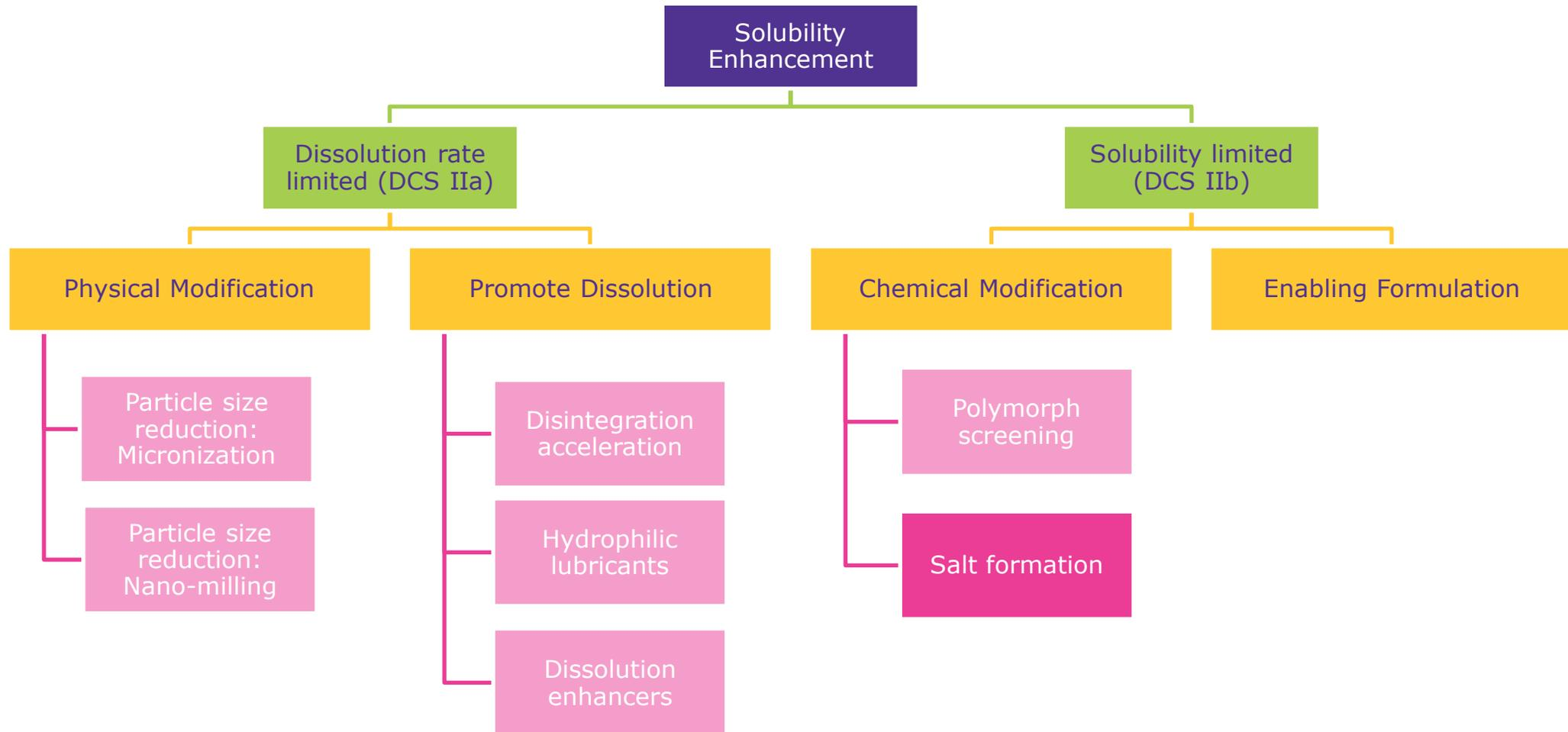


- API polymorphs are a **very important factor for IP protection**:
  - Innovators screen and identify as many polymorphs as possible and protect the related IP, thus **preventing generics entering the market**
  - Non-protected, viable polymorphic forms are **opportunities for generic companies** to enter markets earlier
- Diverse **solvents are broadly used** for polymorph screening, e.g.:

1-Butanol	Chloroform	Ethyl acetate
2-Propanol	Dichloromethane	Formic acid
Acetic acid (glacial)	Diethyl ether	Methanol
Acetone	Dimethyl Sulfoxide	Tetrahydrofuran
Acetonitrile	Ethanol absolute	

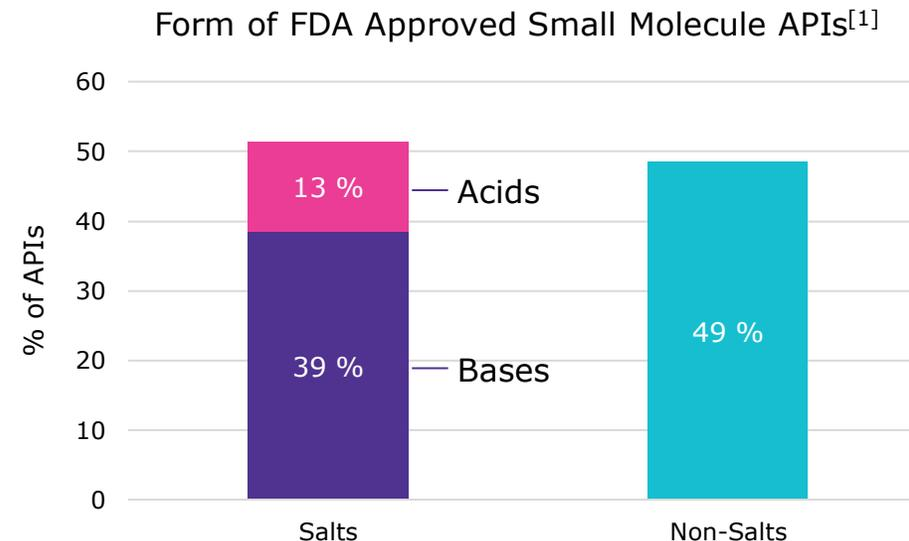
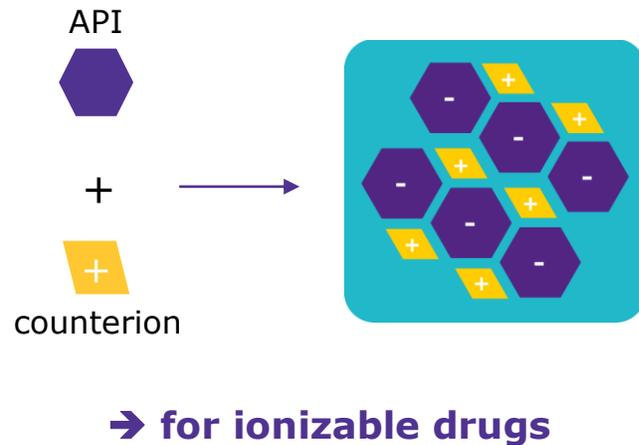


# Strategies and Enabling Technologies for Enhancing API Solubility



# Salt Formation

- Salt formation is a **well-established** technique, employed for several decades
- Currently marketed drugs are **predominantly in salt form** (~50%), still with increasing trend
- **Principle:** API is ionized with the aid of a counterion (acid or base, typically H<sup>+</sup> transfer)



~ 50 %  
of drug molecules  
are administered  
in salt form<sup>[1]</sup>

[1] Trends in Active Pharmaceutical Ingredient Salt Selection based on Analysis of the Orange Book Database, Paulekuhn *et al.*, 2007

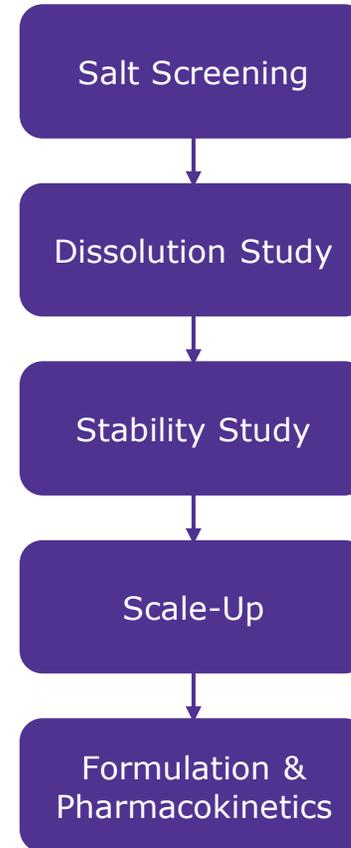


# Salt Formation – Why Salts?

## Opportunity: Optimizing solid state properties

- Enhanced bioavailability<sup>[1]</sup>
  - Solubility
  - Dissolution rate
- May also result in beneficial impact on:
  - API purity
  - Physical & chemical stability
  - Manufacturability
  - Particle size
  - Flowability

## Typical Process Flow



## Benefits of Salts

- Well established
- Simple & cost-effective
- High regulatory acceptance

## Challenges

- API needs to be ionizable
- Common Ion Effect
- Disproportionation
- Hygroscopicity

[1] Use of pharmaceutical salts and cocrystals to address the issue of poor solubility, Holm *et al.*, 2013



# Salt Formation – Examples

## Examples marketed drugs in salt form<sup>[1]</sup>:

Product	API	Counterion	Company
Clindac (antibiotic)	Clindamycin	Phosphate	Sandoz
Baytril (antibiotic)	Enrofloxacin	Mesylate	Bayer
Trimeton (antihistaminic)	Chlorpheniramine	Maleate	MSD
Escitalopram (antidepressant)	Escitalopram	Oxalate	Hexal
Gastrografin (contrast agent)	Diatrizoate	Meglumine	Bayer
Ketorol (anti-inflammatory)	Ketorolac	Thromethamine	Dr. Reddy`s

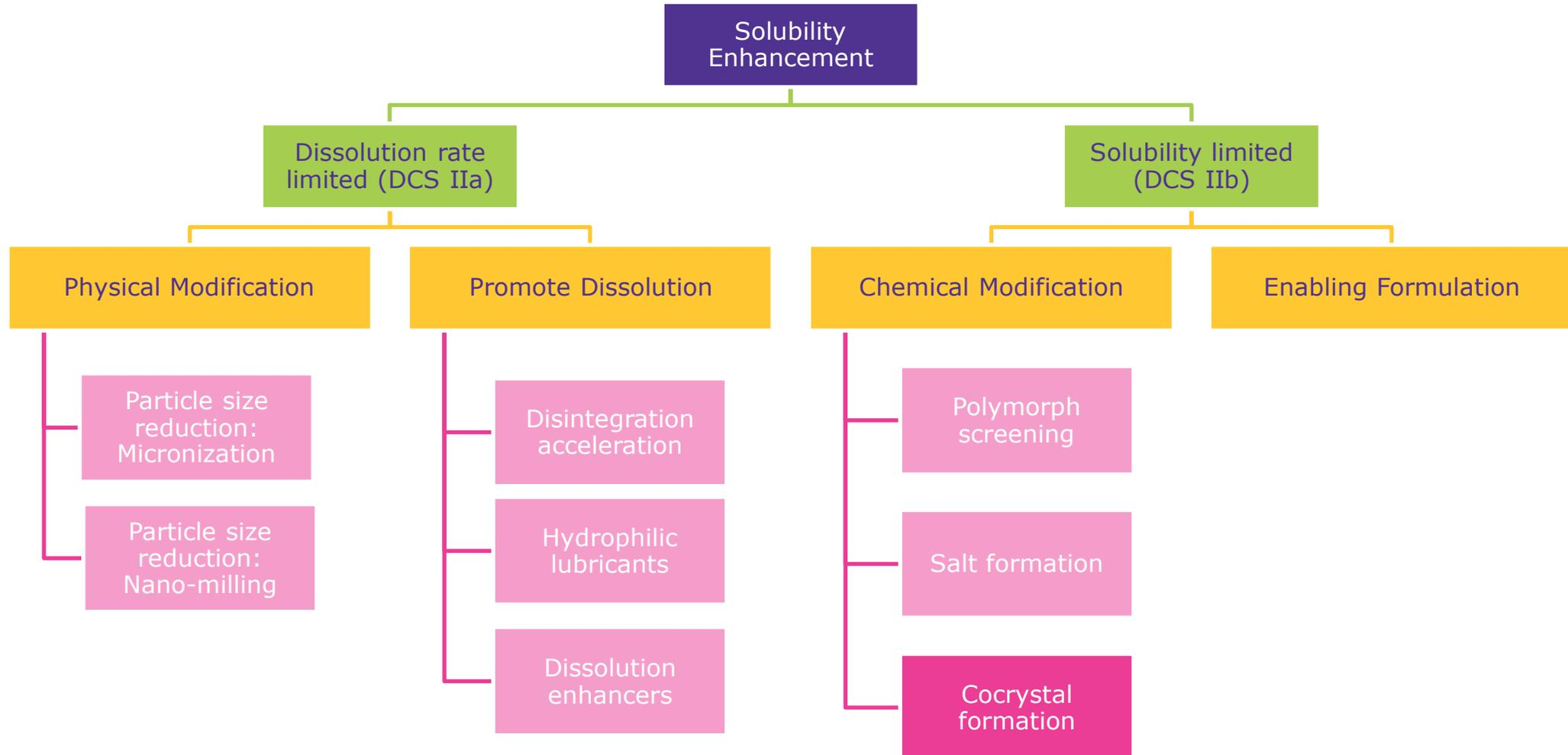
## Common salt former / counterions:

Acetic acid	Citric acid	Imidazole	Salicylic acid
Adipic acid	Formic acid	Lactic acid	Sorbic acid
Benzenesulfonic acid	Fumaric acid	Maleic acid	Succinic acid
Benzoic acid	L-Glutamine	Malic acid	Sulfuric acid
Boric acid	Hydrochloric acid	Ortho-Phosphoric acid	Tartaric acid

[1] Pharmaceutical salts: A summary on doses of salt formers from the Orange Book, Saal & Becker, 2013; [2] Internal Evaluation

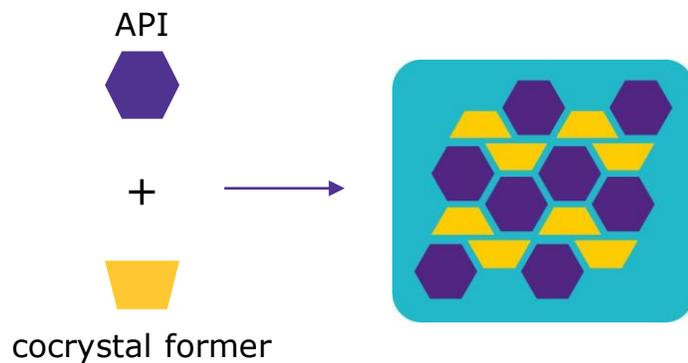


# Strategies and Enabling Technologies for Enhancing API Solubility

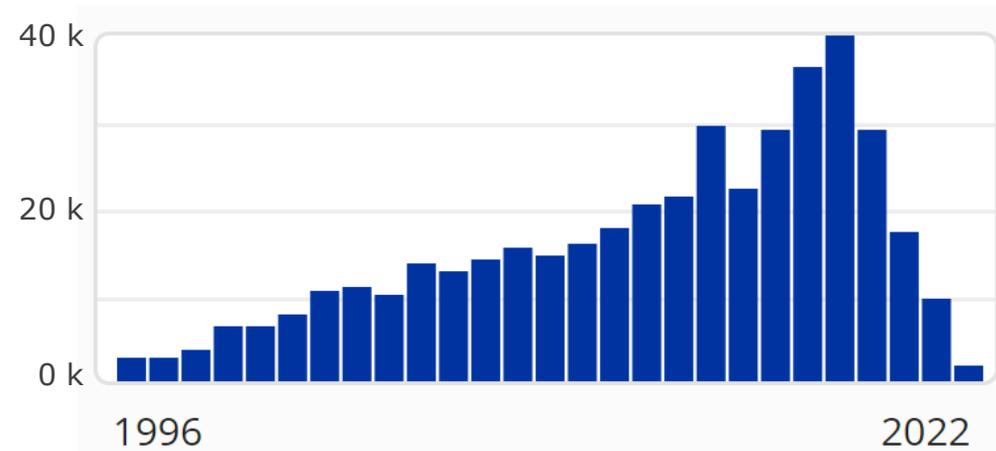


# Cocrystal Formation

- Pharmaceutical cocrystal formation is a **relatively new approach**
- **Alternative for non-ionizable compounds**, as nearly all kinds of APIs can form cocrystals
- **Principle:** API and co-former interact by certain forces and build one joint crystal lattice



→ for non-ionizable drugs



Reference hits for "co-crystals" in Formulus database search (patents applications) per year (2018: 39 k; date 10/22).

Marketed API cocrystals are still rare (<10) but increasing<sup>[1]</sup>

[1] Recent Advances in Pharmaceutical Cocrystals: From Bench to Market, Bandaru *et al.*, 2021

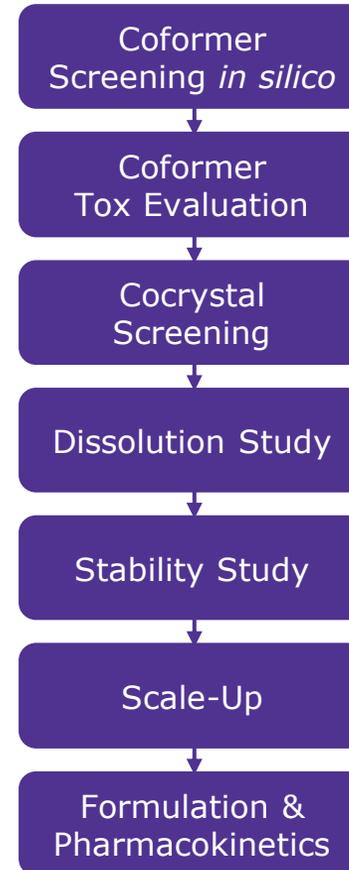


# Cocrystal Formation – Why Cocrystals?

## Opportunity: Optimizing solid state properties

- Enhanced bioavailability<sup>[1,2]</sup>
  - Solubility
  - Dissolution rate
- May also result in beneficial impact on:
  - API purity
  - Physical & chemical stability
  - Manufacturability
  - Particle size
  - Flowability

## Typical Process Flow



## Benefits of Cocrystals

- Alternative if non-ionizable
- No common ion effect
- No disproportionation effects
- Less prone to hygroscopicity

## Challenges

- Coformer selection
- Crystallization process control
- Regulatory acceptance

[1] N. R. Hornedo, S. J. Nehm, A. Jayasankar in Encyclopedia of Pharmaceutical Technology, Third Edition, Informa Healthcare, 2006, pp. 615–635.

[2] N. Upadhyay, T. P. Shukla, A. Mathur, J. S. K. Manmohan, Pharmaceutical Sciences Review and Research 2011, 8, 144–148.



# Cocrystal Formation – Examples

## Examples marketed cocrystal drugs<sup>[1]</sup>:

Product	API : Coformer	Company
Suglat®	Ipragliflozin : L-Proline	Astellas / MSD
Entresto®	Valsartan : Sacubitril	Novartis
Steglatro®	Ertugliflozin : Z-Pyroglutamic acid	MSD
Depakote®	Valproic acid : Valproate sodium	Abbott
Beta chlor®	Chloral hydrate : Betaine	Mead Johnson

## Common cocrystal former:

2,4-dihydroxybenzoic acid	Nicotinamide	L-proline
3,4-dihydroxybenzoic acid	Tert-butyl hydroquinone	Maltol
Tartaric acid	Resorcinol	Oxalic acid
Citric acid	Orotic acid	Urea
Fumaric acid	Sorbic acid	Glycerol
Phloroglucinol	L-tyrosine	Sulfamic acid

## Diverse techniques exist for cocrystal formation:

- Slurrying
- Grinding (solvent-assisted)
- Solvent evaporation
- Sublimation
- Hot melt extrusion (DSC screening)

**Computational co-former screening** is a promising approach to save time and resources

[1] Recent Advances in Pharmaceutical Cocrystals: From Bench to Market, Bandaru et al., 2021; [2] Internal Evaluation



## EMA – Reflection on Cocrystals<sup>[1]</sup>

“[...] There is no strict borderline between the salt formation in the one end with complete proton transfer and cocrystals formation in the other end with no proton transfer. [...] Cocrystals and salts share many conceptual similarities and therefore also similar principles for documentation should be applied.”

## FDA – Guidance on Cocrystals for Industry<sup>[2]</sup>

“[...] Co-crystals are distinguished from salts because unlike salts, the components that co-exist in the co-crystal lattice with a defined stoichiometry interact nonionically. [...] A co-crystal with a pharmaceutically acceptable cofomer [...] can be considered to be a pharmaceutical cocrystal and has a regulatory classification similar to that of a polymorph of the API. Specifically, it is not regarded as a new API.”

## Regulatory Implications of Cocrystals

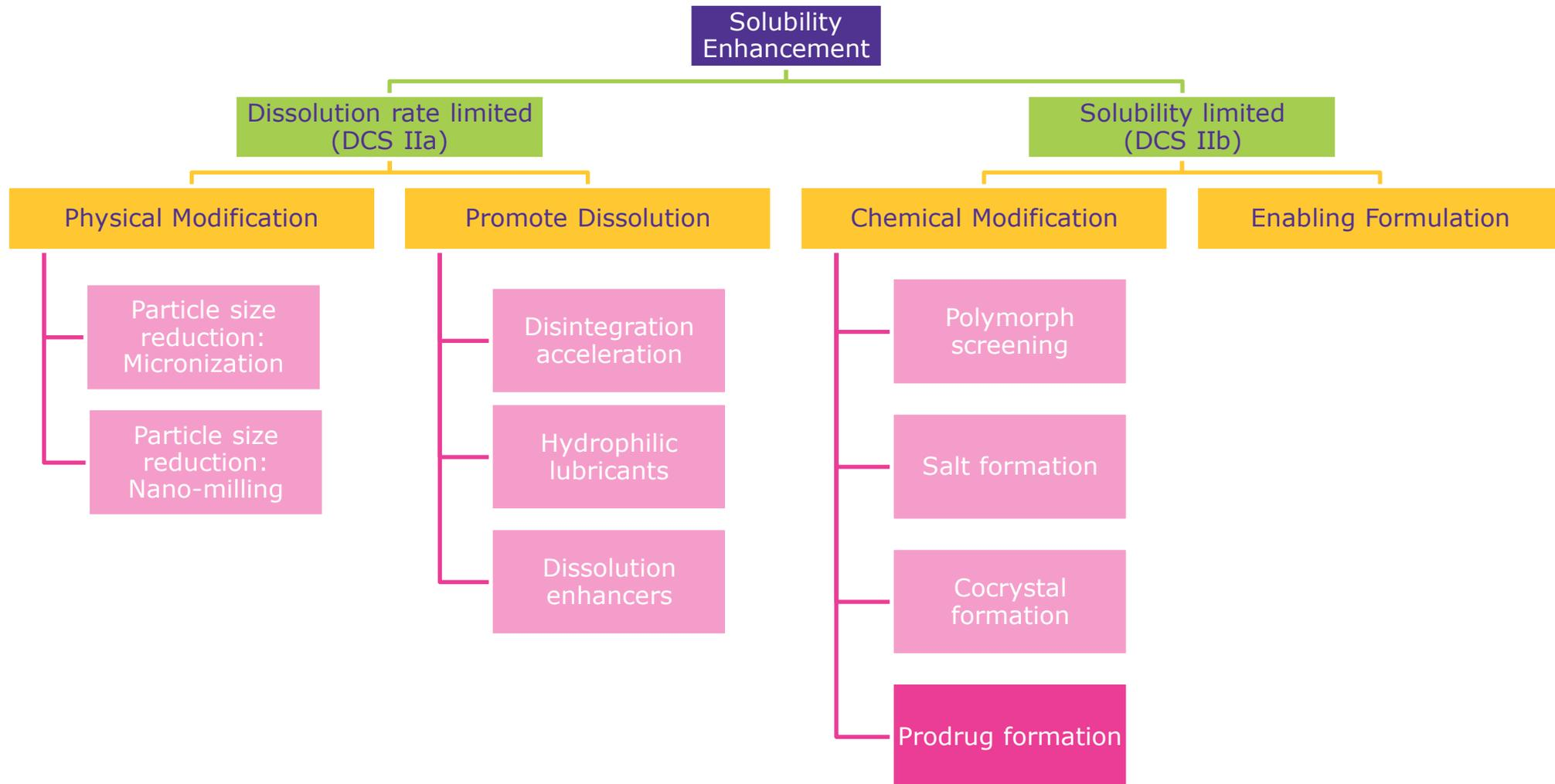
- Reflections on cocrystals (still) strongly differ between FDA and EMA
- Both FDA and EMA provide little guidance regarding co-former

[1] Reflection paper on the use of cocrystals of active substances in medicinal products, EMA, 2015

[2] Regulatory Classification of Pharmaceutical Co-Crystals – Guidance for Industry, FDA, 2018



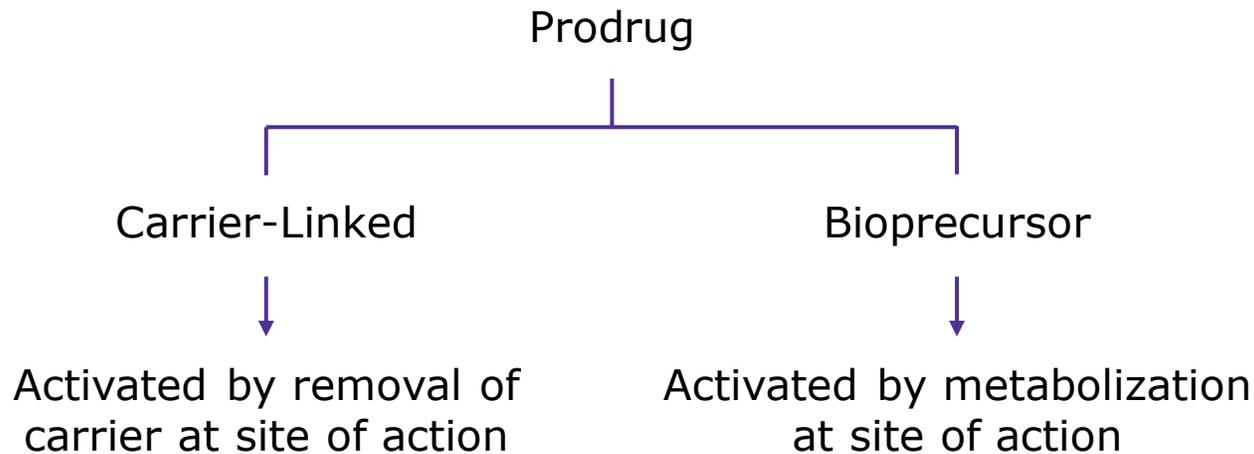
# Strategies and Enabling Technologies for Enhancing API Solubility



# Prodrug Formation

- **Prodrugs** are molecularly modified APIs with improved physicochemical characteristics
- They have no or **low activity until converted** to the active parent drug *in vivo*
- Often regarded as “strategy of last resort”, but offers **various opportunities**<sup>[1]</sup>

Approx. 10 % of all marketed drugs are prodrugs – with increasing trend<sup>[2]</sup>



## Objectives of prodrug approach:

- Improved solubility
- Targeted release
- Improved membrane permeability
- Reduced side effects

[1] The Prodrug Approach: A Successful Tool for Improving Drug Solubility, FDA, 2015

[2] The expanding role of prodrugs in contemporary drug design and development, Rautio *et al.*, 2018



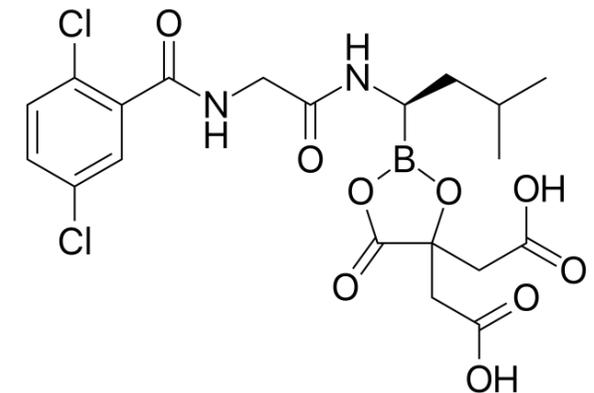
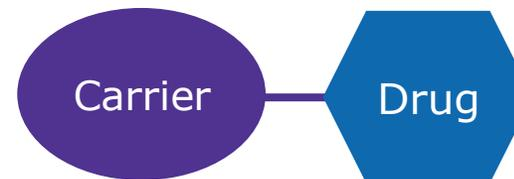
# Prodrug Formation

## Carrier-Linked Prodrugs

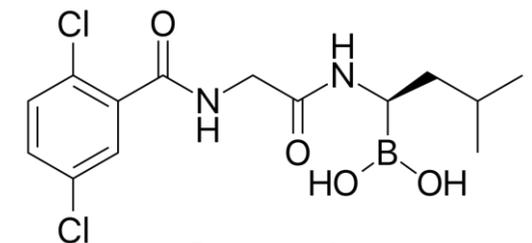
- Common carriers are e.g.: Polymers (e.g. PEGs), proteins, antibodies
- Linkage of carrier to drug possible by diverse covalent binding types, e.g.: ester, amide, carbamate, carbonate, ether, imine, phosphate etc.

### Example: Ixazomib (Takeda, approval 2015<sup>[2]</sup>)

- Indication: Multiple myeloma, reversible proteasome inhibitor
- Prodrug hydrolyzes under physiological conditions to active form



Ixazomib Citrate (Prodrug)



Ixazomib

[1] Prodrug approach: An overview of recent cases, Abet *et al.*, 2017

[2] The expanding role of prodrugs in contemporary drug design and development, Rautio *et al.*, 2018

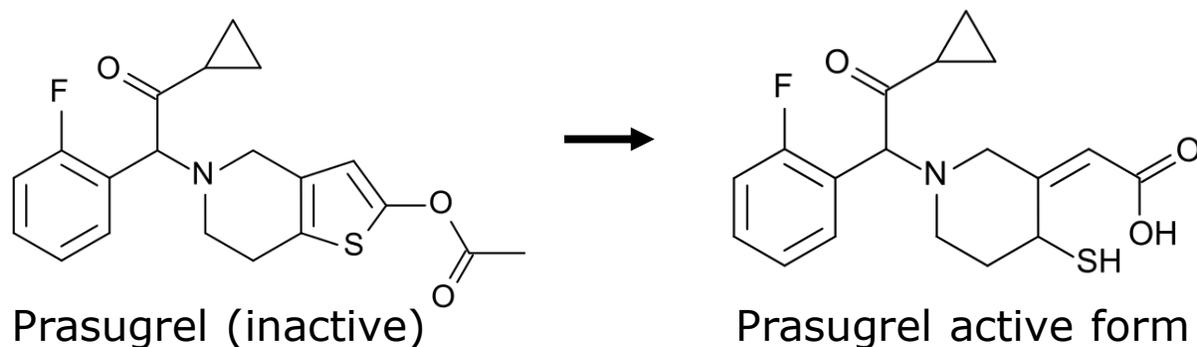


## Bioprecursors

- No carrier or linker, precursor acts as substrate to enzymatic activation<sup>[1]</sup>
- Activation by diverse chemical modification possible, e.g. reductive or oxidative activation, phosphorylation, decarboxylation, etc.

**Example:** Prasugrel (Daiichi Sankyo, Eli Lilly, approval 2009<sup>[2]</sup>)

- Indication: Thrombotic and cardiovascular events



## Benefits of Prodrugs

- Versatile approach
- Can improve several properties
- Alternative if other strategies are not feasible

## Challenges

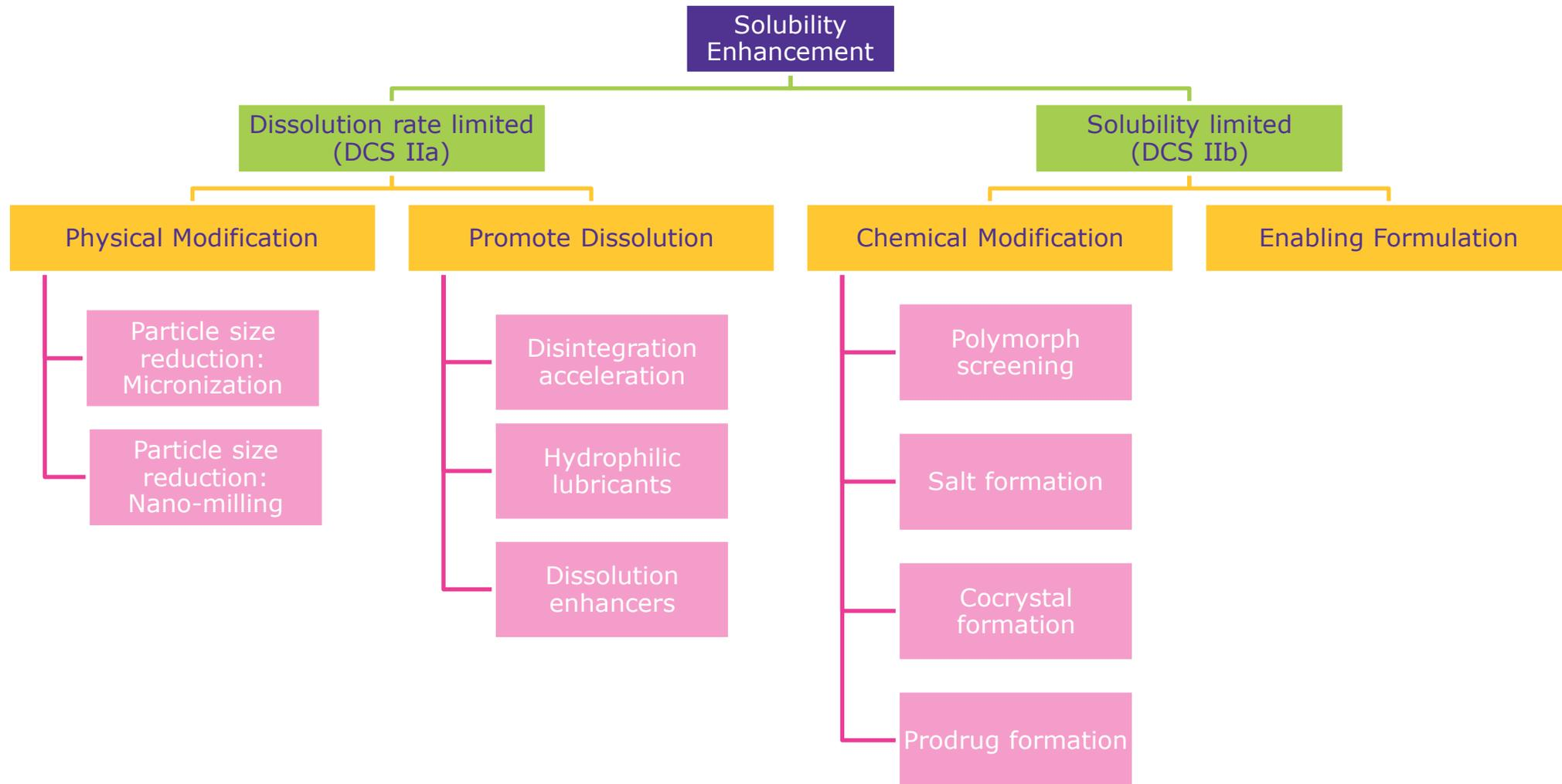
- Demanding development
- Cleaved groups must be safe

[1] Prodrug approach: An overview of recent cases, Abet *et al.*, 2017

[2] Pharmacokinetics and pharmacodynamics of prasugrel, a thienopyridine P2Y<sub>12</sub> inhibitor, Dobesh, 2009

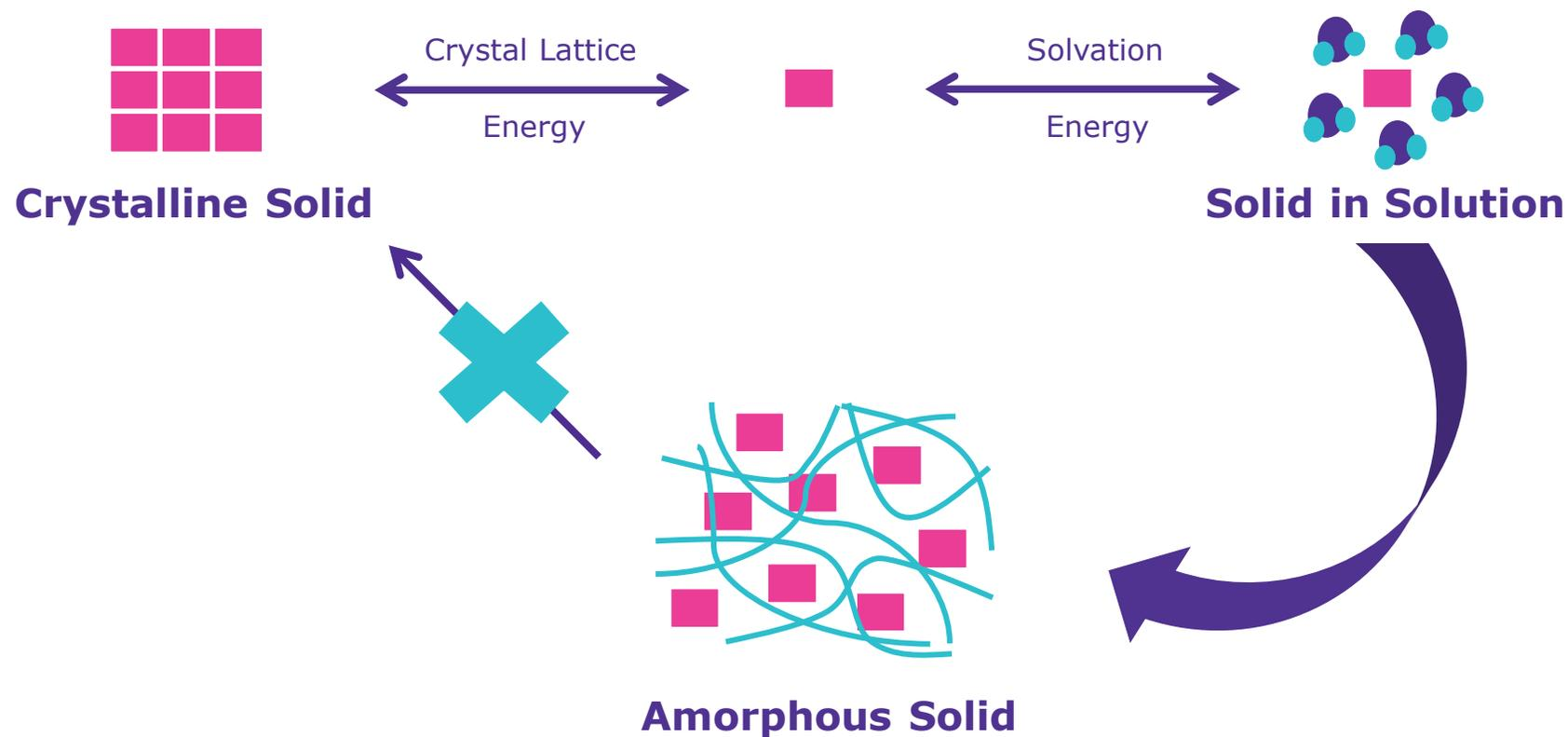


# Strategies and Enabling Technologies for Enhancing API Solubility



# Solid dispersion technologies - Solid state modification

## Solubility of DCS IIb molecules can be enhanced via the amorphous form

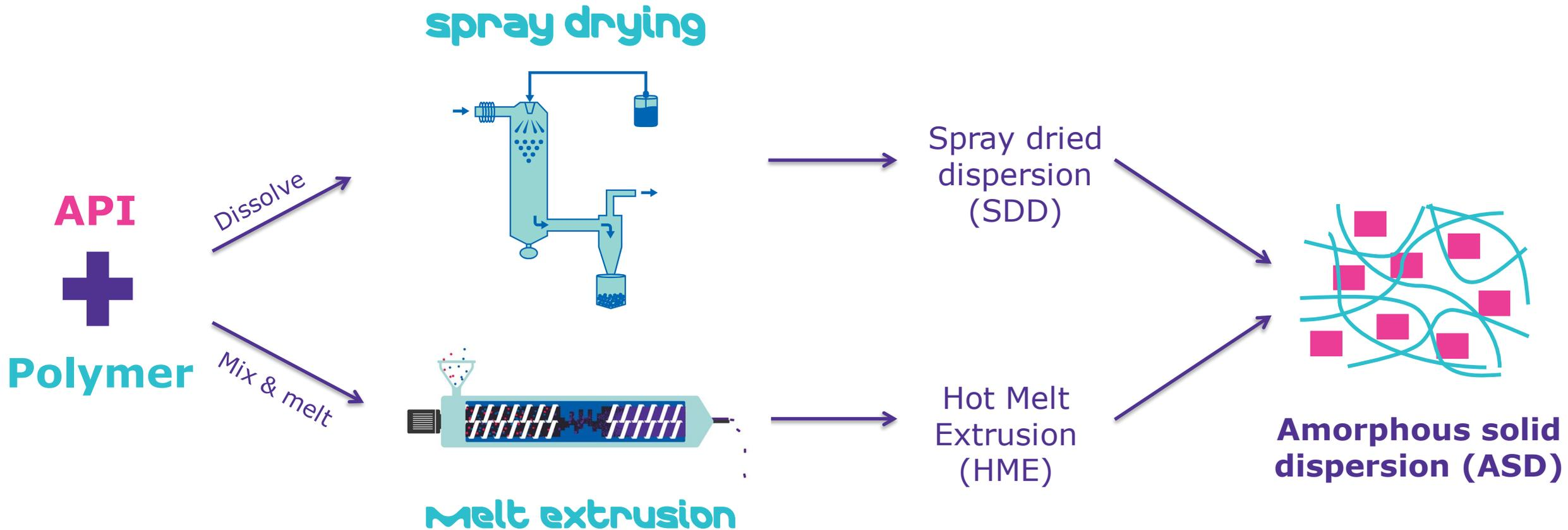


**>> The amorphous solid-state can be leveraged to enhance solubility of DCS IIb**



# Amorphous solid dispersions

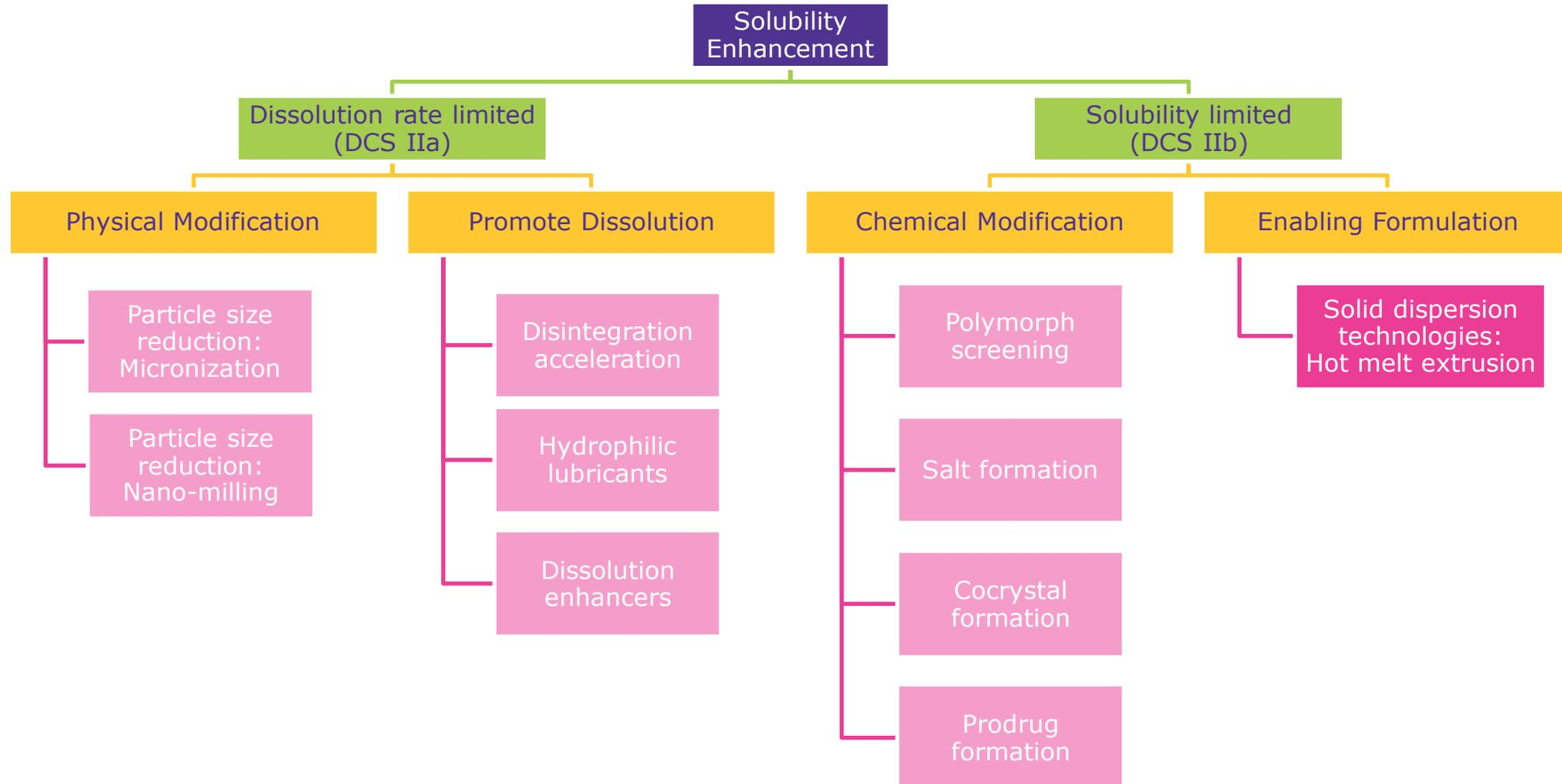
The most common amorphous formulation have a polymeric matrix



Goal is a homogenous dispersion of drug and polymer



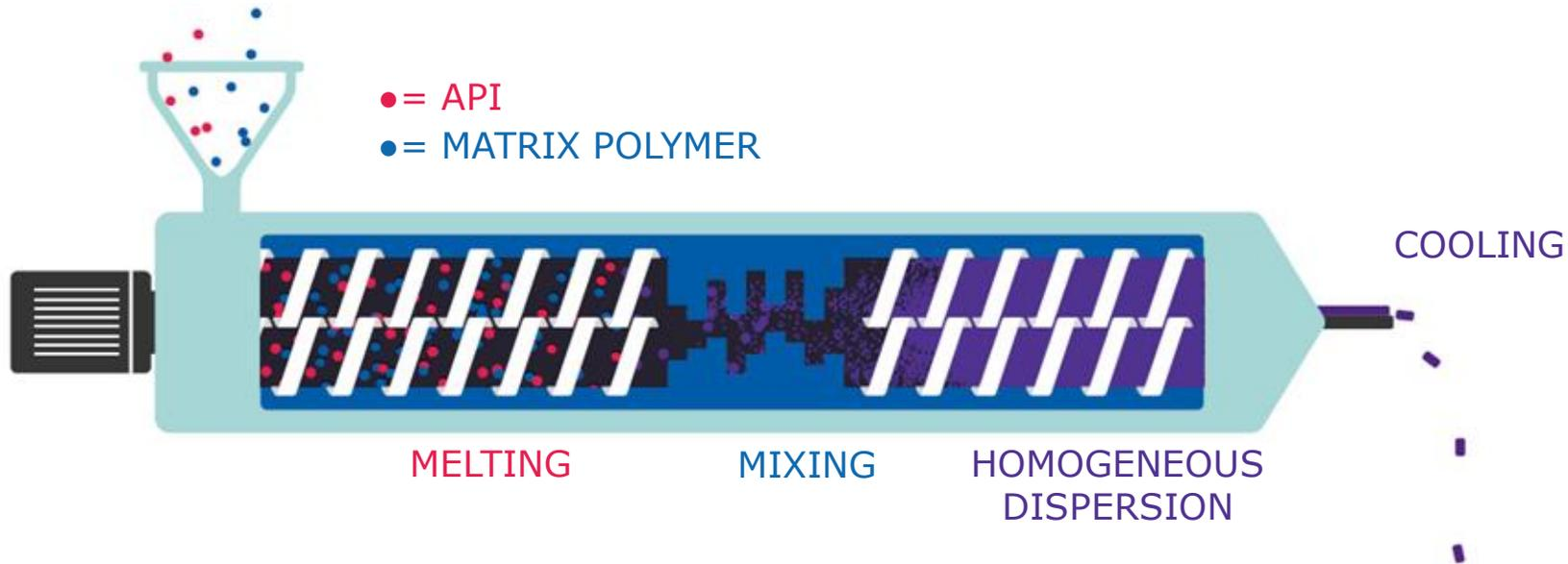
# Strategies and Enabling Technologies for Enhancing API Solubility



# Amorphous solid dispersions

## Advantages of the HME process

SAFC®



### Advantages of HME

- Enhanced solubility & bioavailability
- Stable, high drug loading
- Controlled release
- Continuous effective process
- Solvent-free manufacturing
- Support of various dosage forms

 **MIX. MELT. PERFORM.**

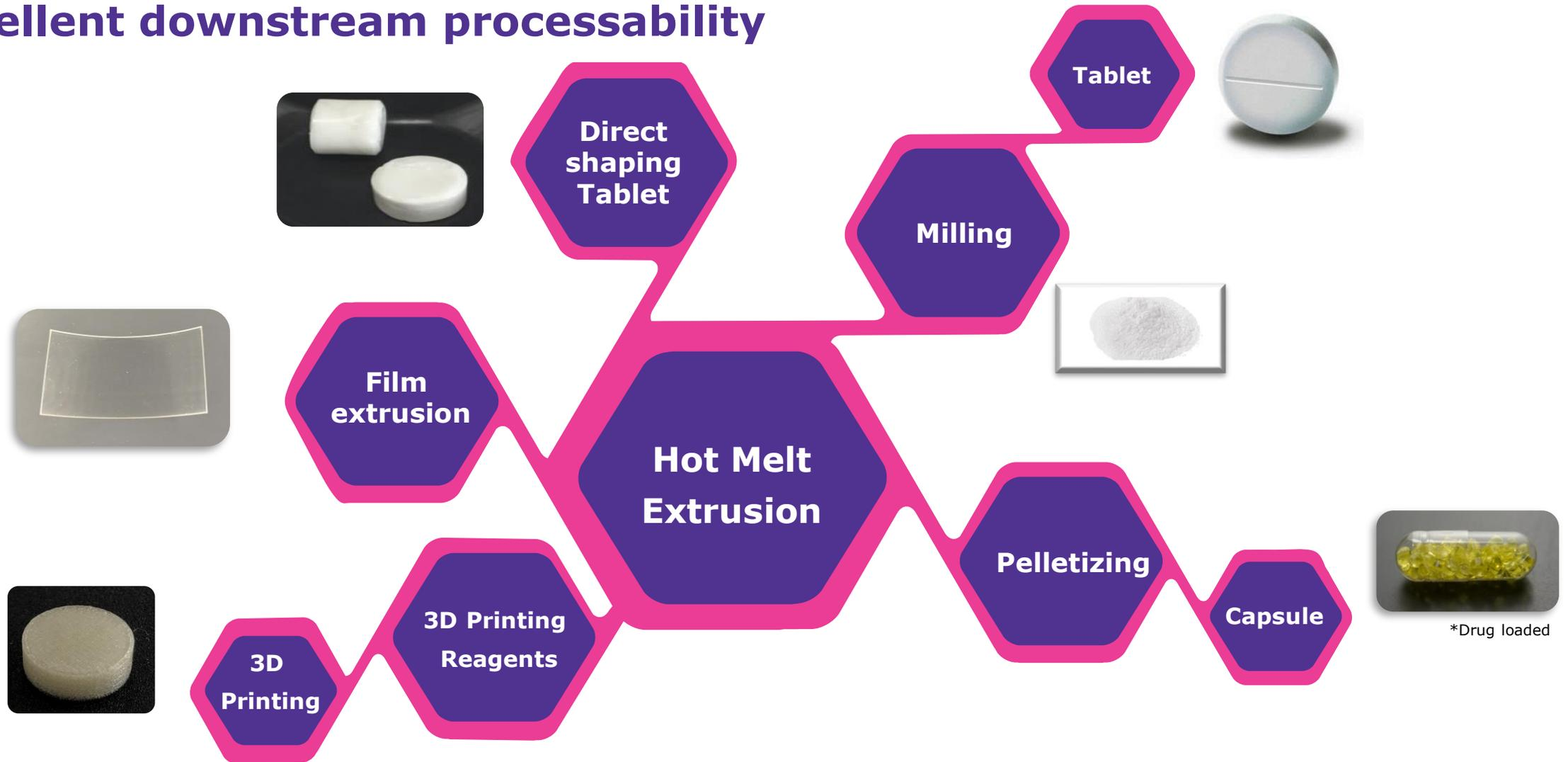




# PVA for Hot Melt Extrusion – Downstream flexibility

## Excellent downstream processability

SAFC®



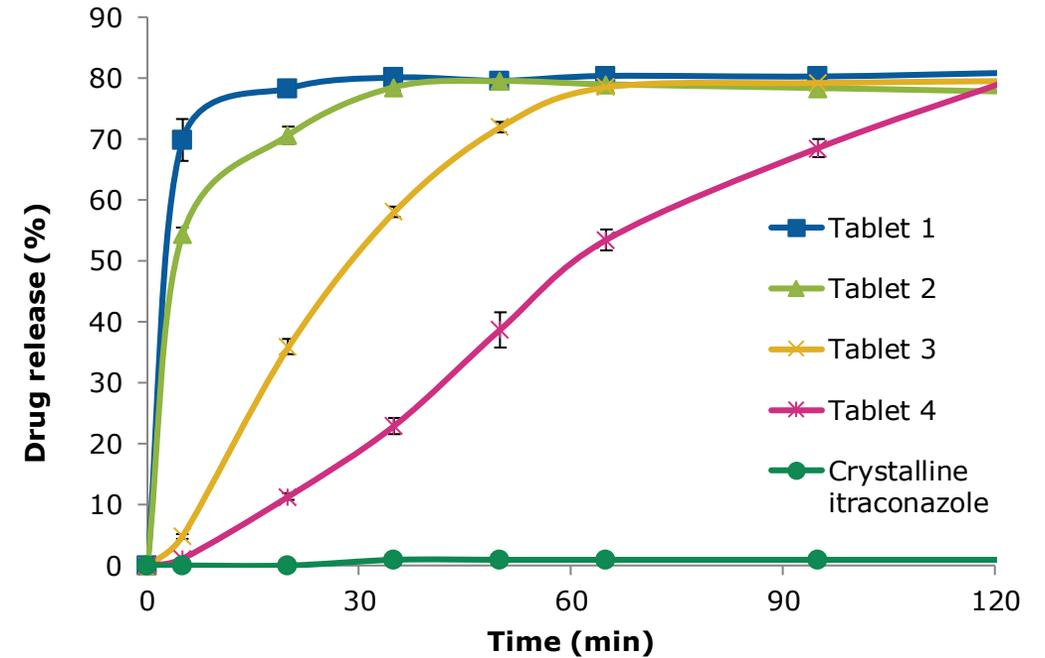
**Versatile polymer extrudates can be processed into various dosage forms**



# PVA for Hot Melt Extrusion – Release profile

## Addition of other excipients allow flexible release profiles

	Tablet 1	Tablet 2	Tablet 3	Tablet 4
Extrudate (%)	50	50	50	60
Microcrystalline cellulose (%)	10	10	10	10
K <sub>2</sub> CO <sub>3</sub> (%)	-	-	14.75	10
NaCl (%)	14.75	14.75	-	-
Magnesium stearate (%)	0.5	0.5	0.5	0.5
Lactose (%)	16.25	16.25	16.25	11
Silica (%)	1	1	1	1
Crospovidone (%)	7.5	7.5	7.5	7.5
Compressed force (kN)	15	10	10	10
T <sub>max</sub> (min)	15	30	60	120



Dissolution of compressed tablets based on milled itraconazole:PVA (FDA-recommended conditions for itraconazole, 900 mL SGF, 37 °C, 100 rpm, formulated extrudate with 30 % drug load; n=3)



Single polymer, single extrudate, many options to modify dissolution kinetics

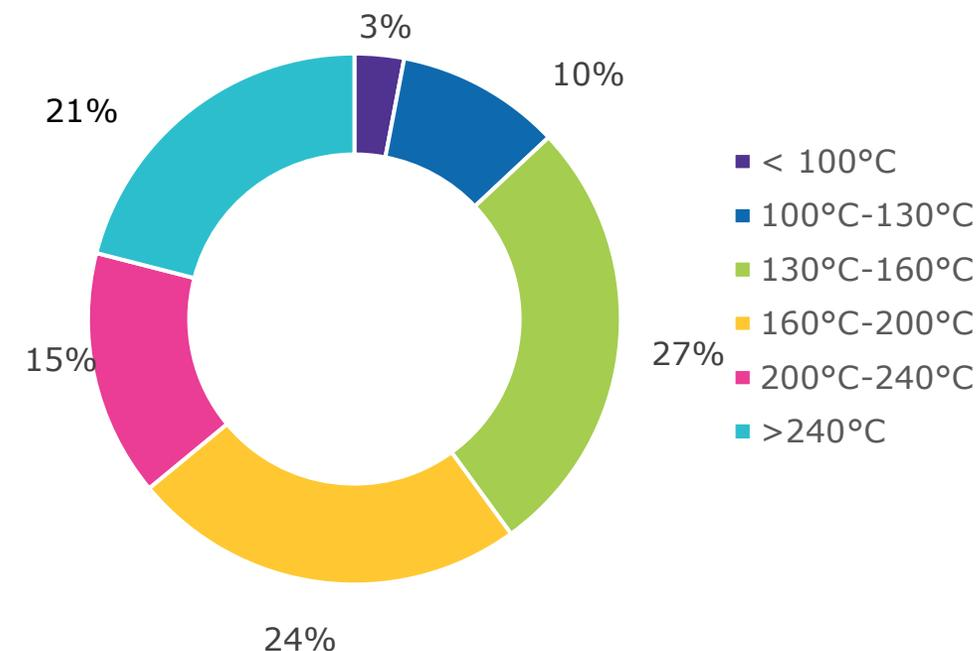


# PVA for Hot Melt Extrusion – Hydrolysis optimization (PVA 4-88)

## Parateck® MXP 4-88: Thermo-stabile polymer for high melting point APIs

API BCS II&IV	T <sub>m</sub> of API	Loading Capacity	Solubility Enhancement (max.)
Ibuprofen*	78 °C	> 30 %	2 x
Cinnarizine	118-122 °C	< 20 %	10 x
Indomethacin	151 °C	> 50 %	3 x
Ketoconazole	146 °C	> 35 %	17 x
Naproxen	152 °C	> 30 %	4 x
Atorvastatin	159-160 °C	> 55 %	154 x
Itraconazole	167 °C	> 30 %	80 x
Carbamazepine	204 °C	> 30 %	2 x
Telmisartan*	260 °C	> 15 %	35 x

\*Plasticizer required



T<sub>m</sub> Breakdown of 67 BCS II and IV compounds:  
 Sarah Shugarts and Leslie Z. Benet, The Role of Transporters in the Pharmacokinetics of Orally Administered Drugs, Expert Review, Pharmaceutical Research, Vol. 26, No. 9, September 2009



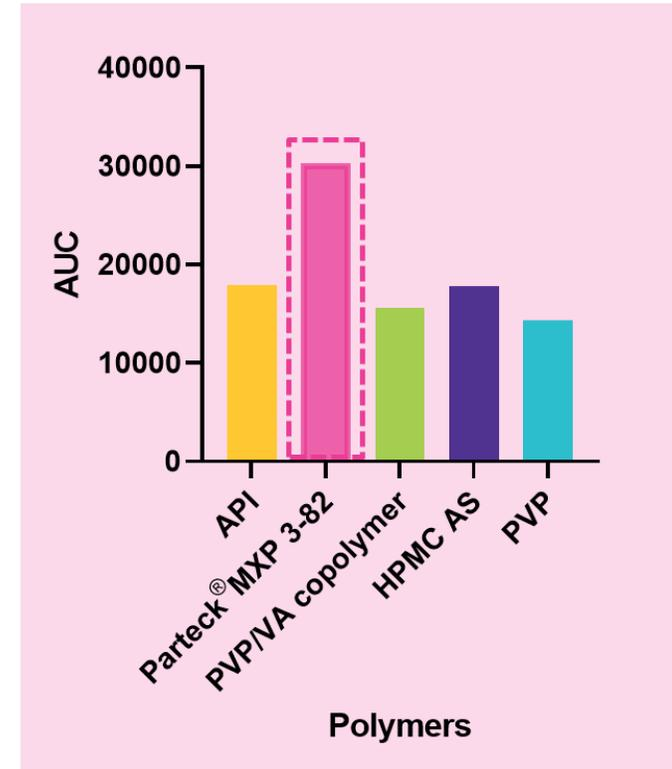
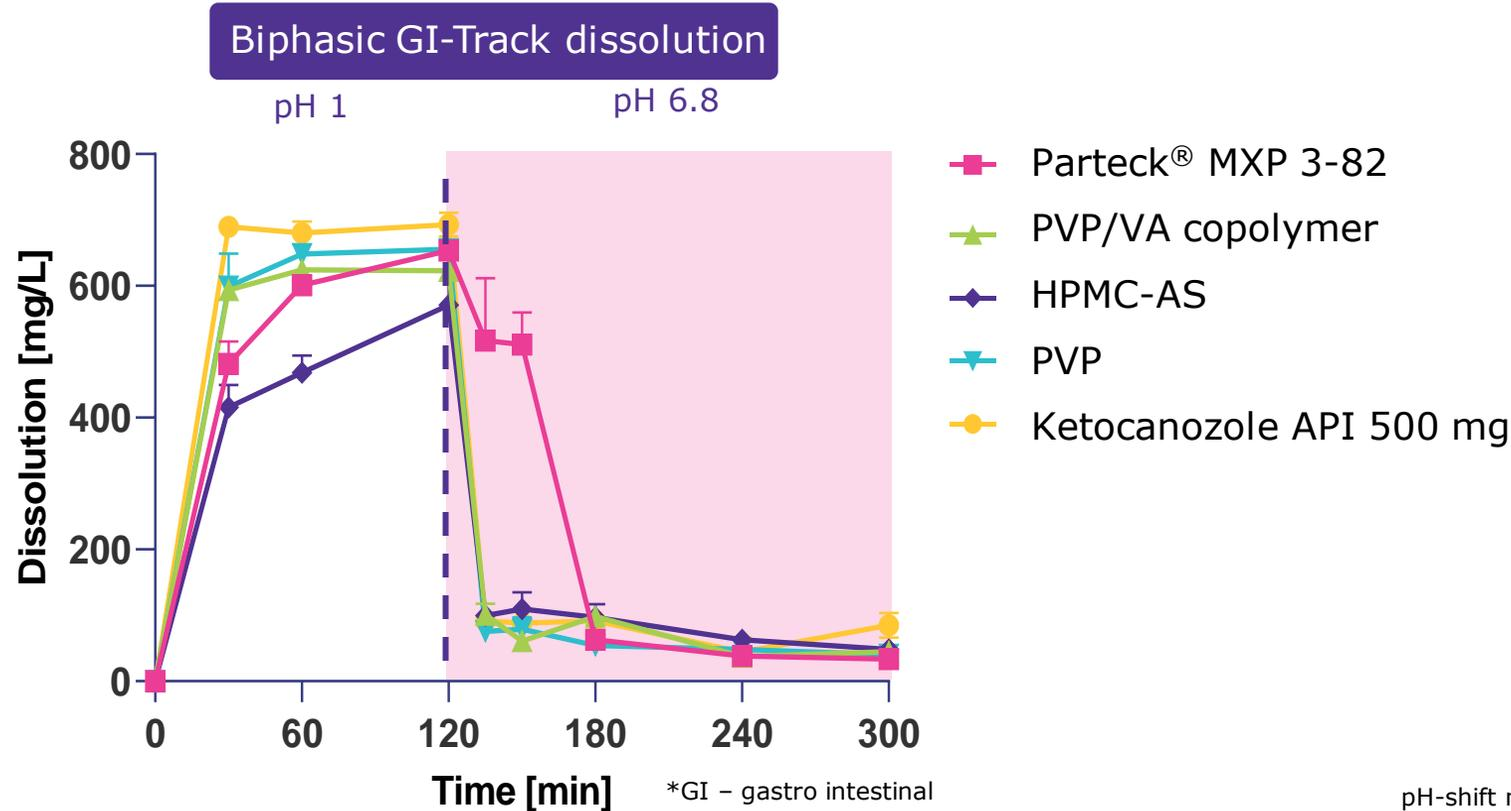
APIs have a range of melting points:

**Good thermo-stability** is needed to cover a wide range (PVA 4-88)



# PVA for Hot Melt Extrusion – Hydrolysis optimization (PVA 3-82)

## Par-teck® MXP 3-82: Formulation performance, prolonged supersaturation of Ketoconazole



pH-shift method (750ml 0.1M HCl for 120 min, add 250ml PP pH 6.8 final 1000ml) HME formulation with 20% Ketoconazole, Paddle: 50 rpm, n=3

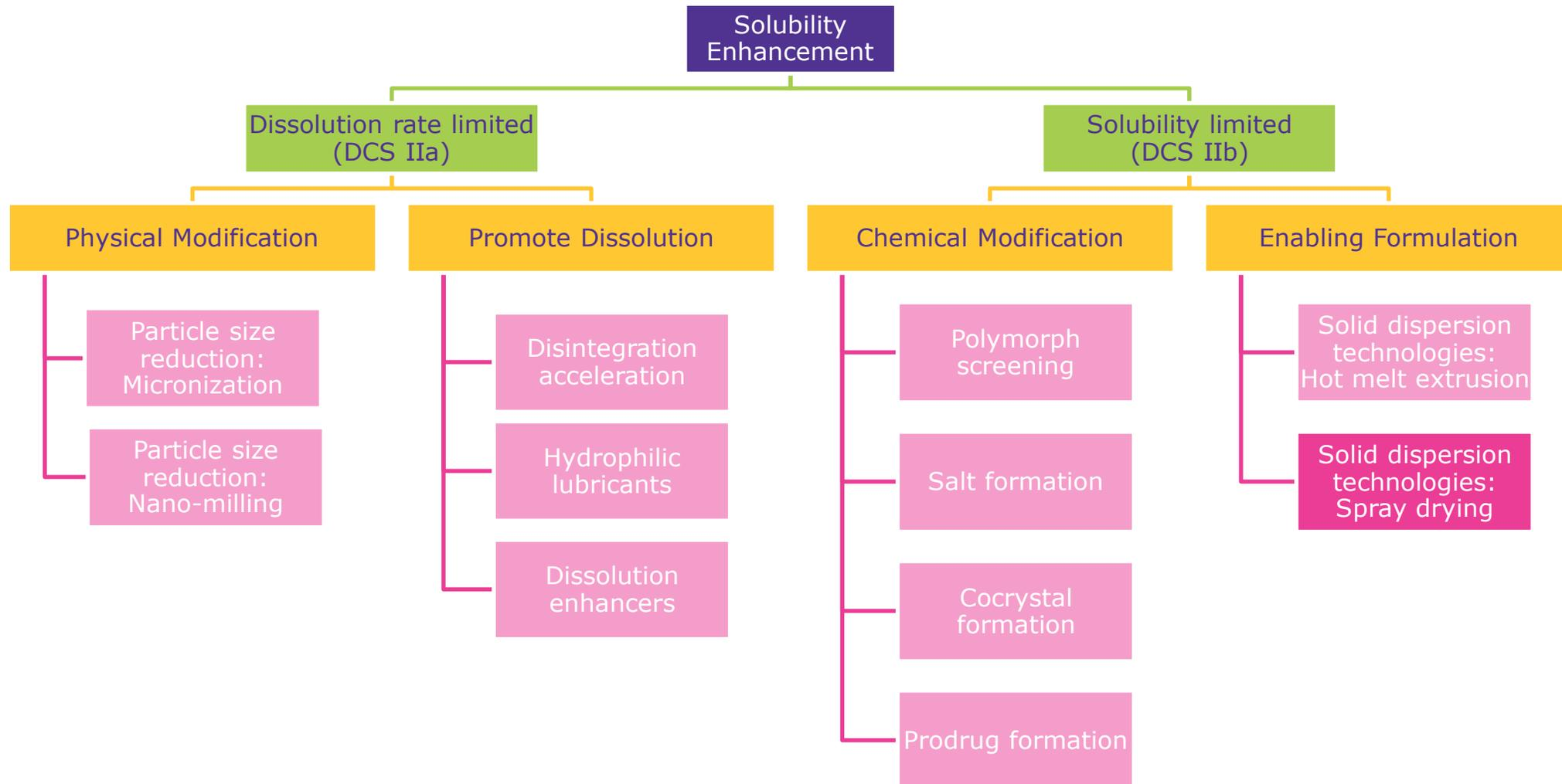
\*AUC – area under the curve



Improved hydrophobicity leads to **superior supersaturation maintenance** via **precipitation inhibition** and prolongates the release profile (PVA 3-82)

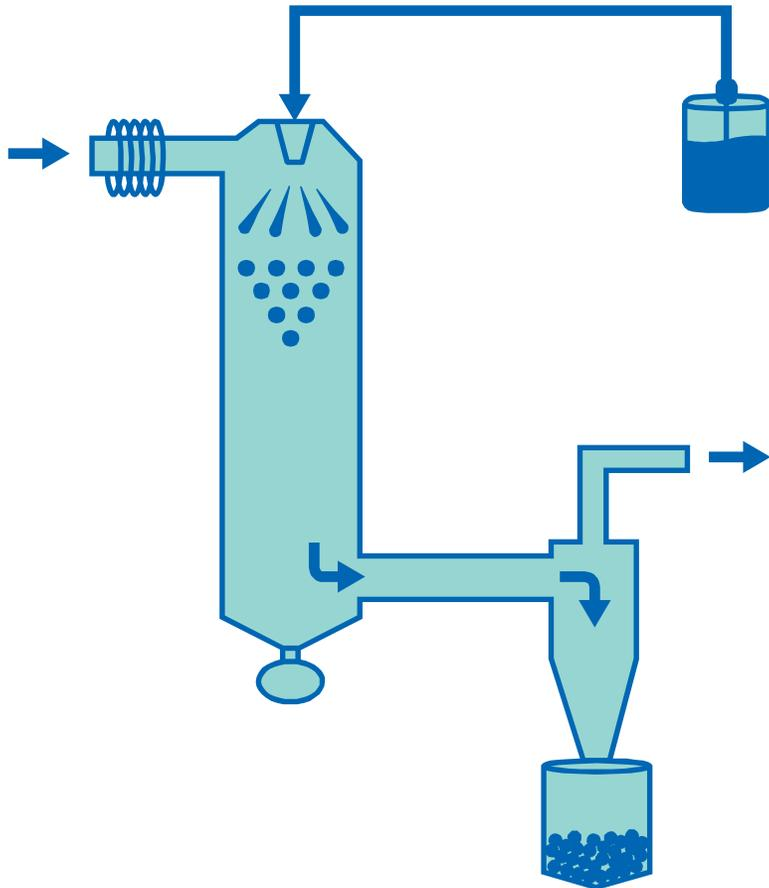


# Strategies and Enabling Technologies for Enhancing API Solubility



# Amorphous solid dispersions

## Advantages of spray drying



### Advantages of spray drying

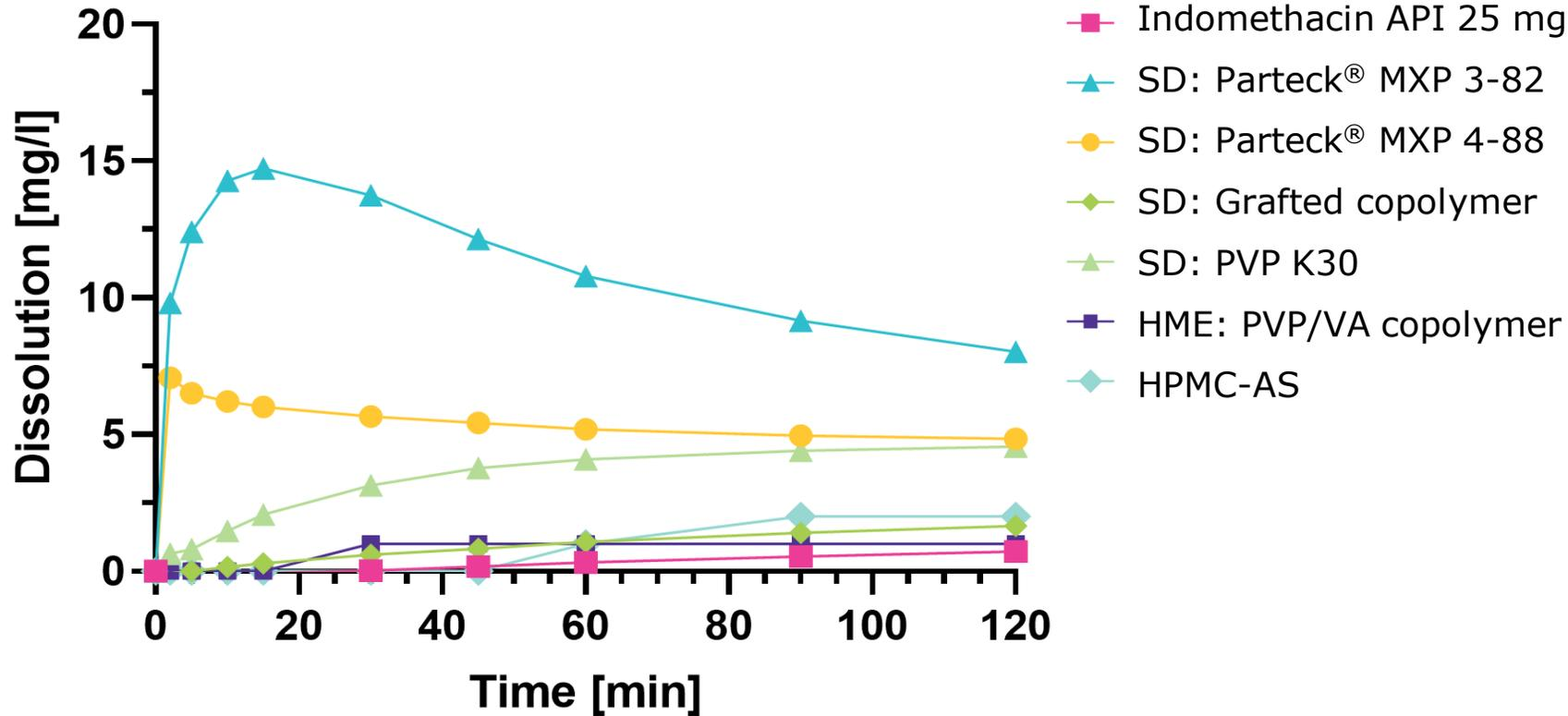
- Enhanced solubility & bioavailability
- Low API amount needed
- Heat sensitive APIs
- Upscaling & high throughput
- Continuous technologies
- Particle engineering

**dissolve. spray. PERFORM.**



# PVA for spray drying

## Dissolution of Indomethacin with an ASD of PVA



- ✓ Fast spring
- ✓ Extended parachute
- ✓ Superior performance

### SD process, Büchi B295

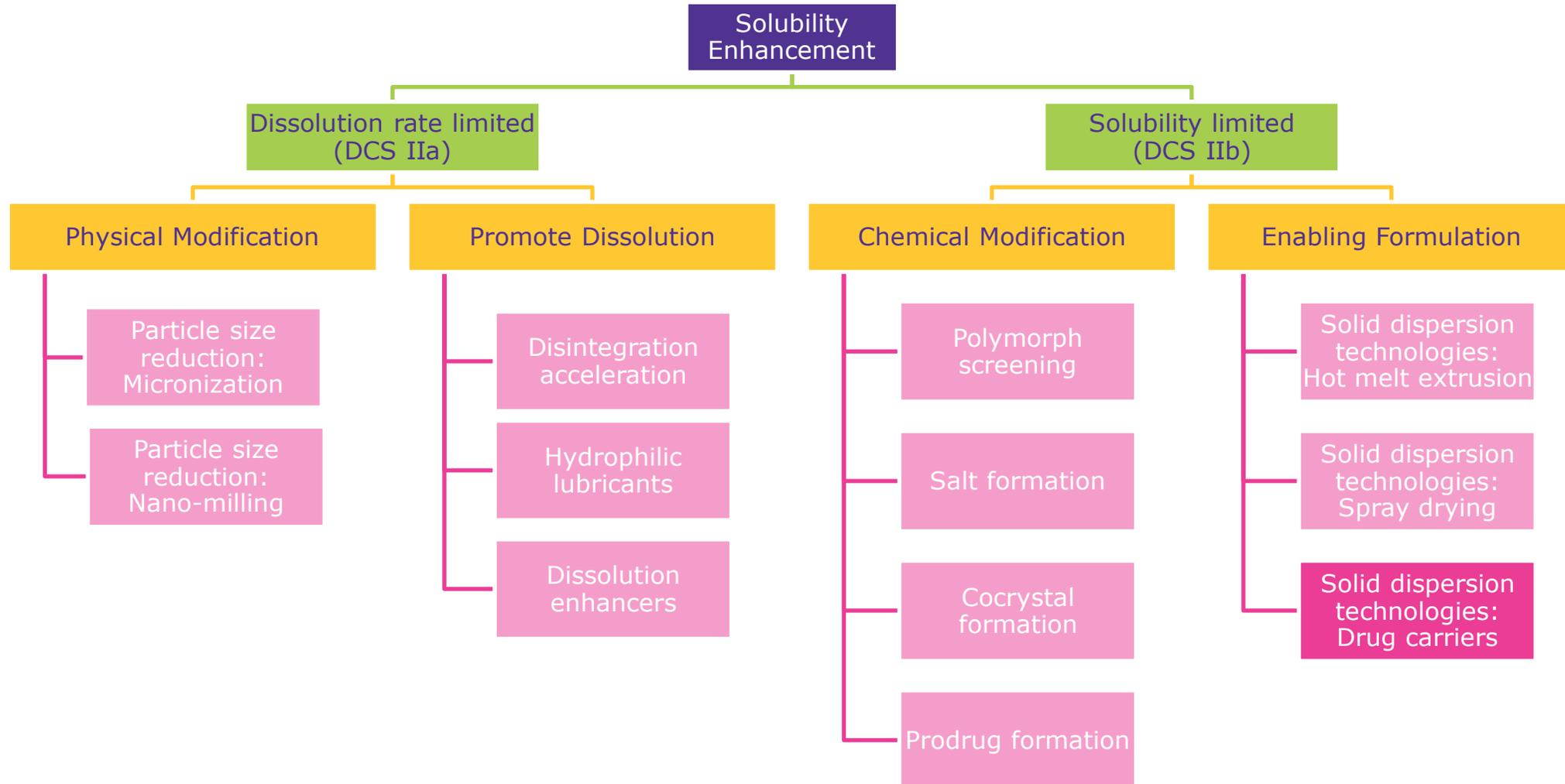
SD via two/three-fluid nozzle: Inlet temperature of 90°C, Outlet temperature of 50°C  
Drying air (N<sub>2</sub>) flow rate of 35 m<sup>3</sup>/h (Aspirator: 100%) Atomization air flow rate of 670 L/min (N<sub>2</sub>: 55mm) Dissolution: Indometacine SGF with 30% DL



**Excellent dissolution enhancement & performance**



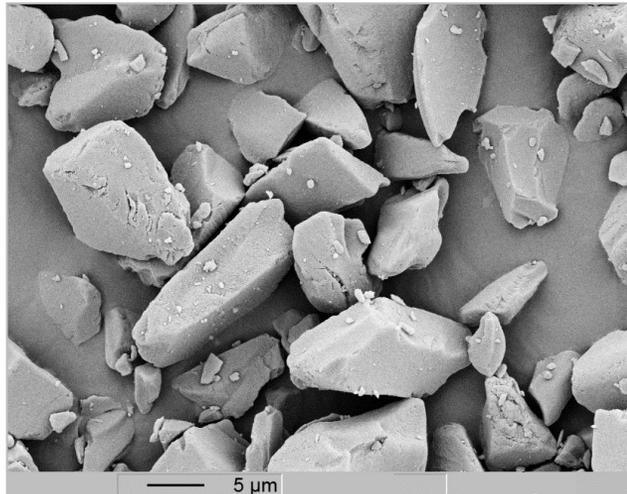
# Strategies and Enabling Technologies for Enhancing API Solubility



# Carrier for amorphous formulation – Mesoporous silica

## Stabilize the amorphous solid-state within the pores of silica

- **Chemical formula:** SiO<sub>2</sub>
- **Pharmacopoeial monograph:** Silicon Dioxide (USP) and Silica, colloidal hydrated (Ph Eur)
- **Regulatory status:** Generally Regarded As Safe (GRAS)\*

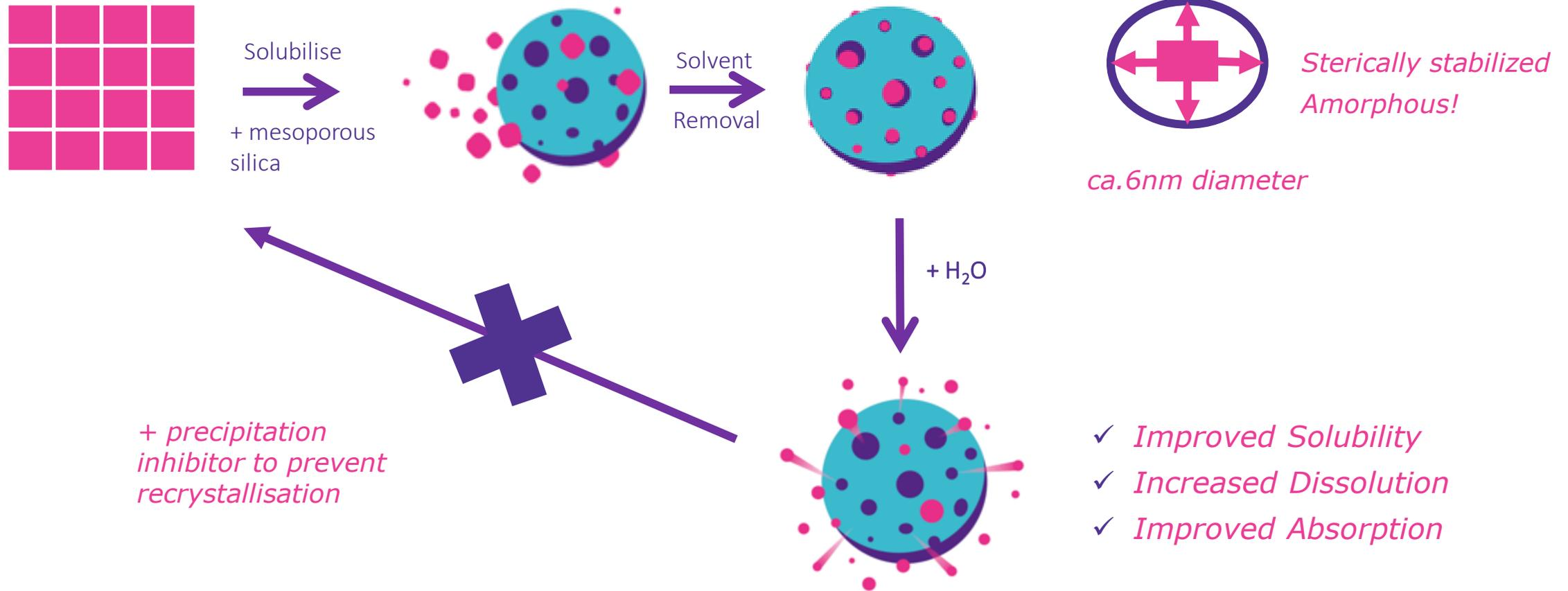


Typical values	
Particle size	5 – 20 μm
Bulk density	0.32 g/mL (0.56 g/mL**)
Surface area	~ 500 m <sup>2</sup> /g
Pore size	~ 6 nm (disordered)



# Mesoporous silica for amorphous formulation

## Stabilize the amorphous solid-state within the pores of silica



 **Pore adsorption and nanoconfinement are the key steps for formulation with mesoporous silica**



# Mesoporous silica for amorphous formulation

## Lab-scale loading of APIs onto mesoporous silica

### Impregnation method



Overhead stirrer

Canulla (API solution)

Vacuum nitrogen

Silica powder

### Suspension method



+ Silica powder  
↓  
API solution

⇒ Stirring  
⇒ Evaporation

 **NO** specialized equipment

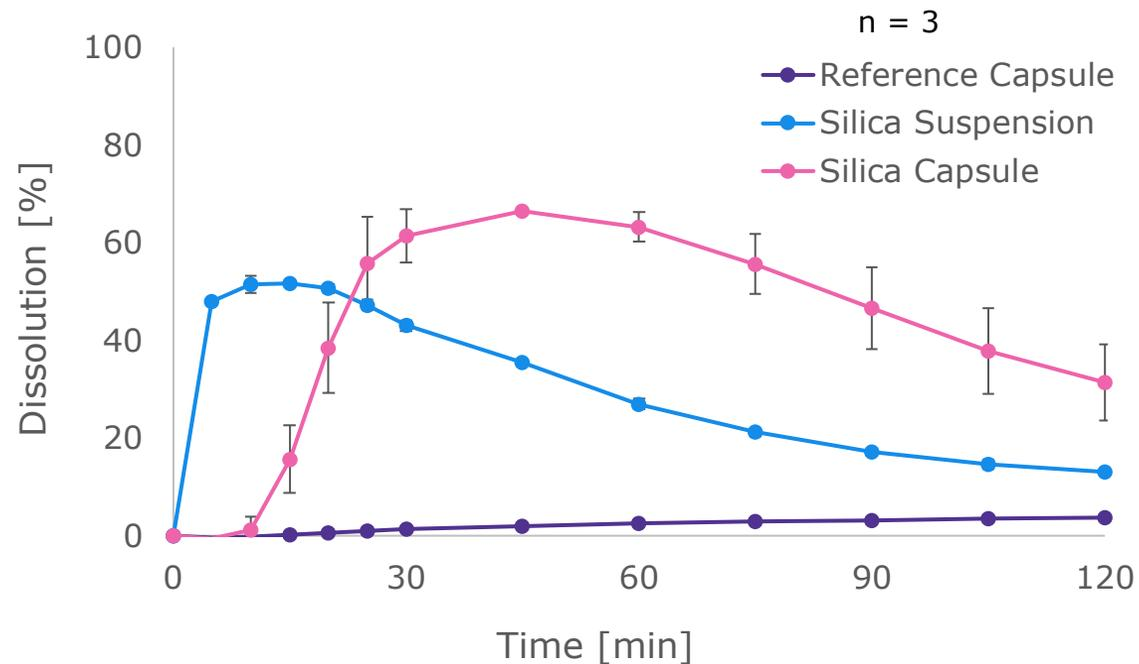
 **NO** extra capital investment



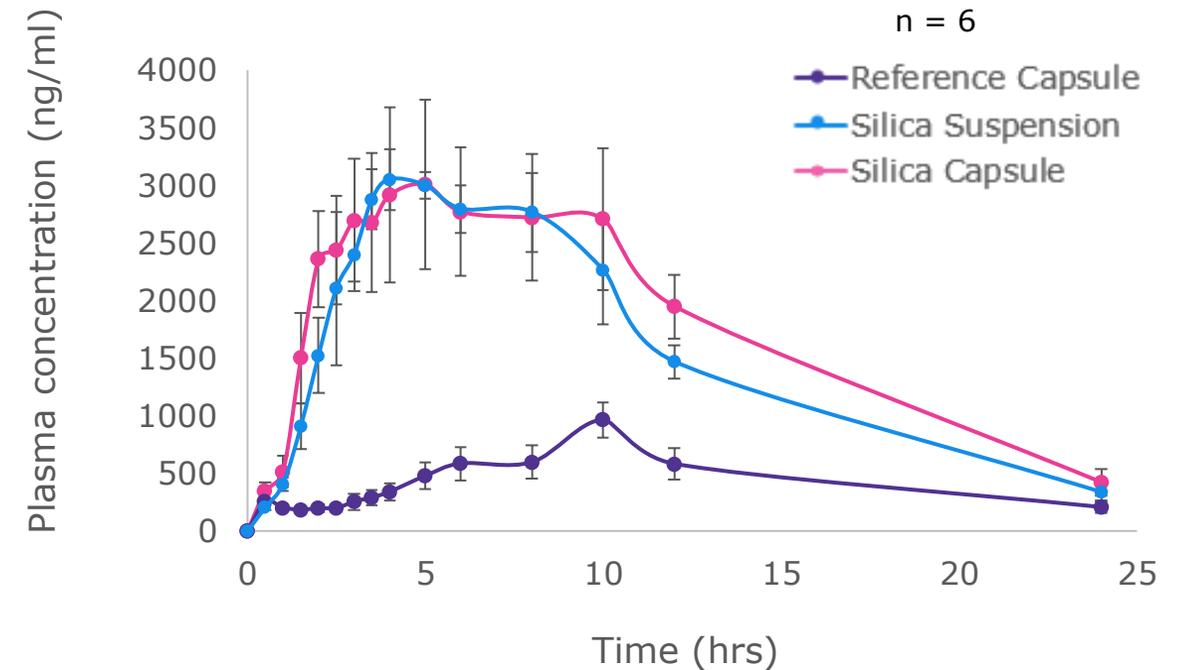
# Mesoporous silica for amorphous formulation

## *In-vitro* and *in-vivo* dissolution profile of fenofibrate

### Biorelevant *in-vitro* dissolution



### *In-vivo* bioavailability in fasted pigs



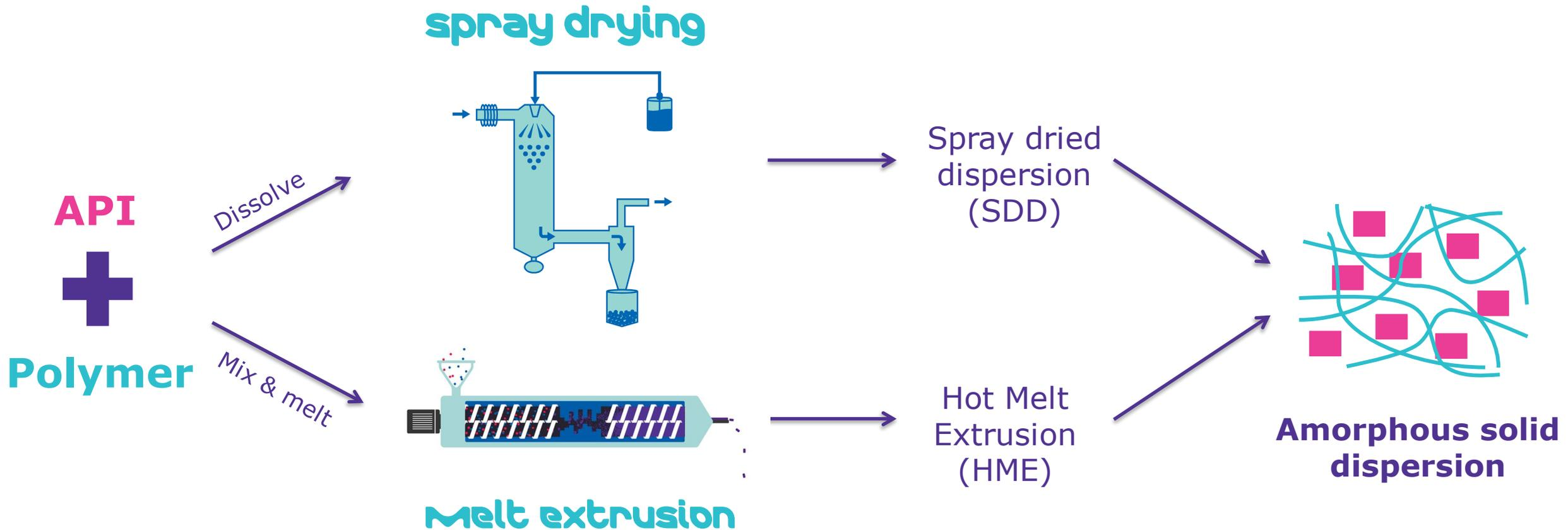
 **Biorelevant dissolution provides a good prediction of relative bioavailability**  
**PK study in pigs indicates a significant bioavailability enhancement**



Amorphous stability

SAFC®

# Polymeric ASDs can lead to phase separation and recrystallization



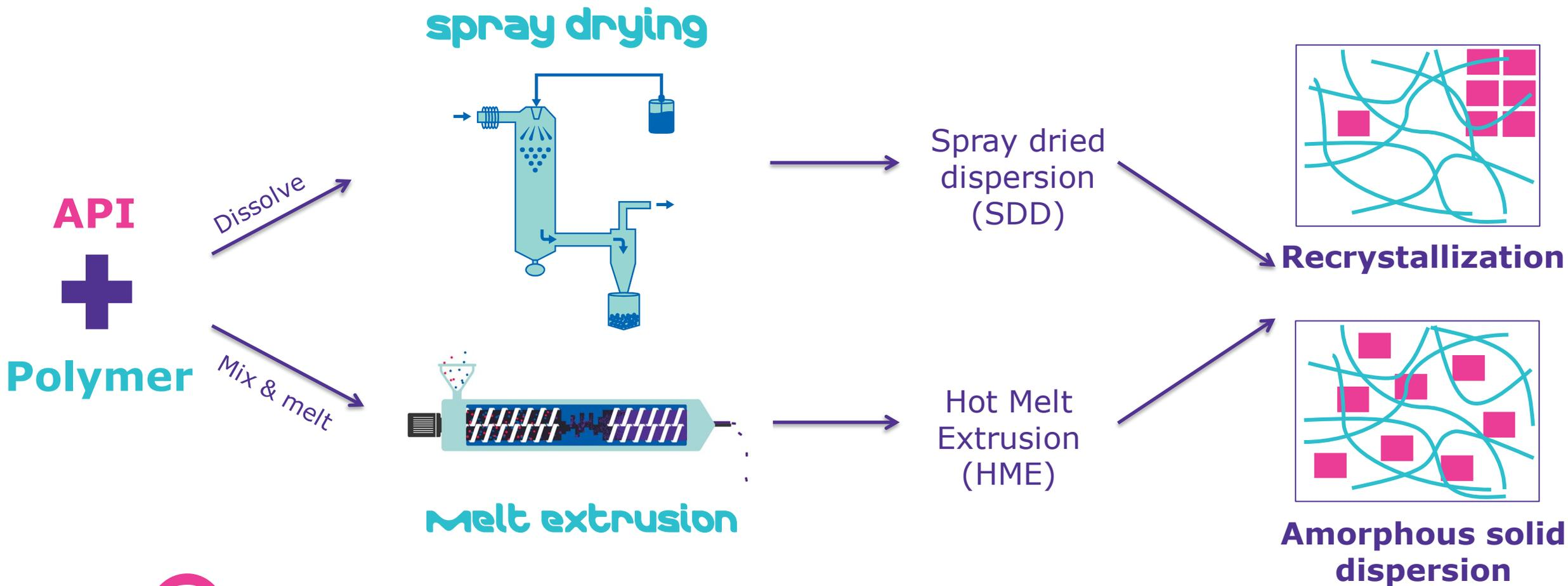
**Goal is a homogenous dispersion of drug and polymer**



Amorphous stability

# Polymeric ASDs can lead to phase separation and recrystallization

SAFC®



How can we predict if our drug will be unstable?

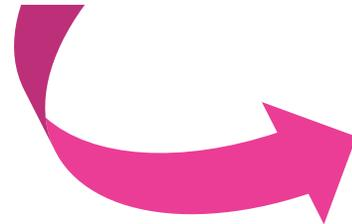


Amorphous stability

## The Glass Forming Ability (GFA) describes the stability in the amorphous form



Poor glass formers (GFA-I) have a higher propensity for re-crystallization. They are more fragile in the amorphous form.



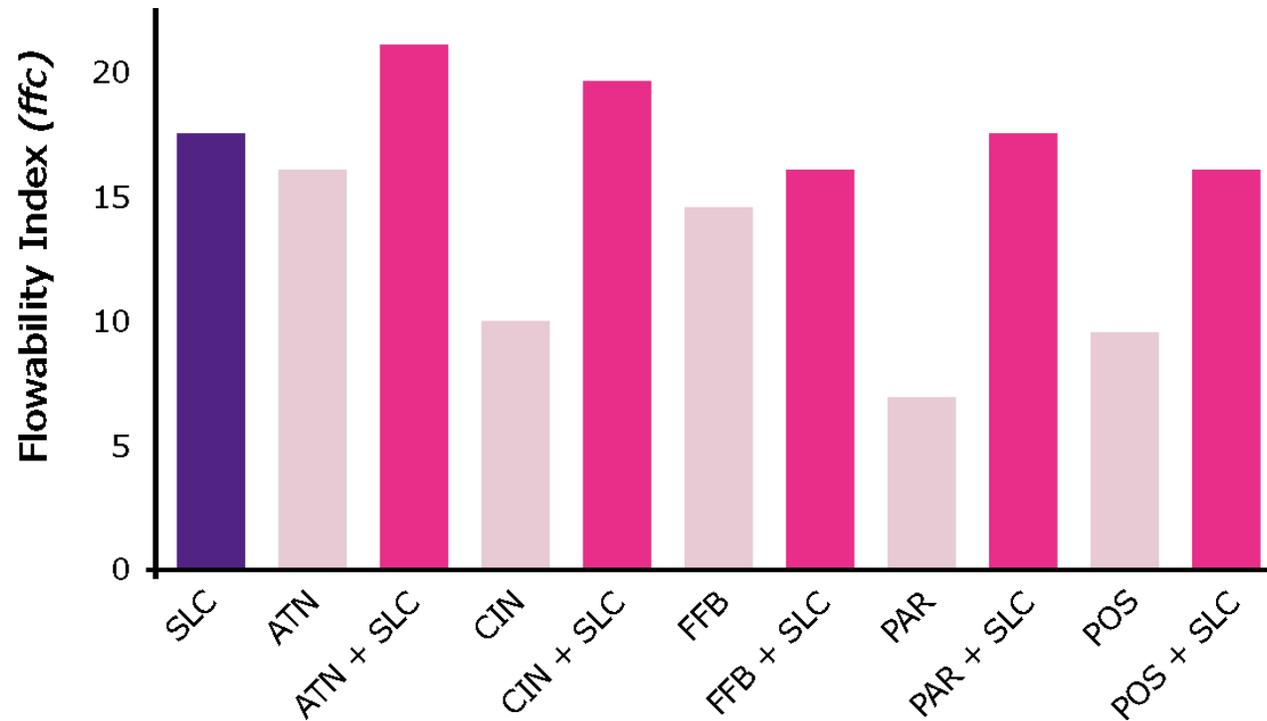
**Mesoporous silica is the best-in-class technology to stabilize poor glass formers**

\*



# Mesoporous silica for amorphous formulation

## Platform technology to stabilize unstable amorphous APIs



### Platform technology approach

#### Uniformity & consistency:

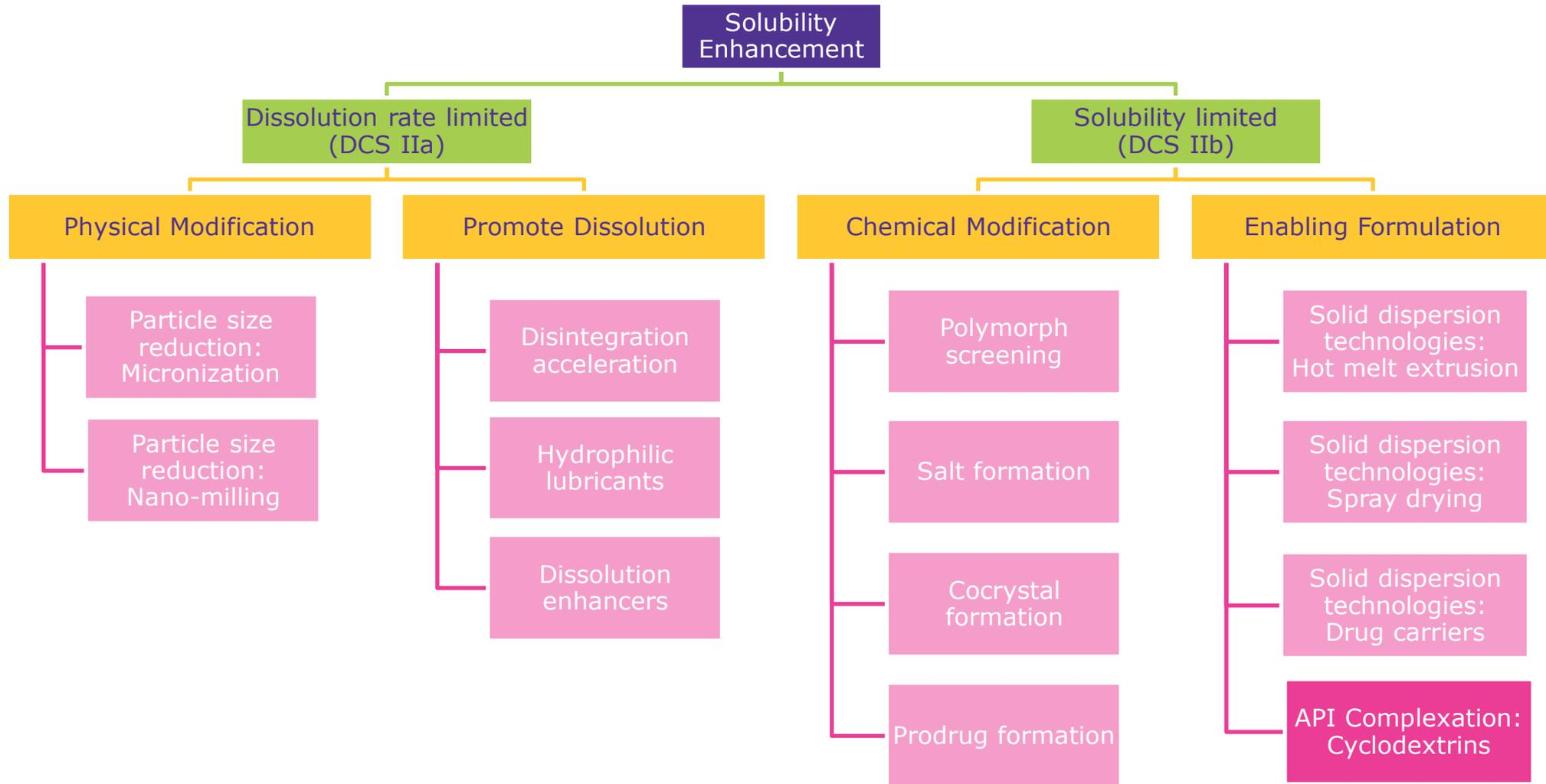
- Particle size
- Flowability
- Tablet hardness
- Friability



De-risk unexpected polymorphic changes and challenges in drug development



# Strategies and Enabling Technologies for Enhancing API Solubility

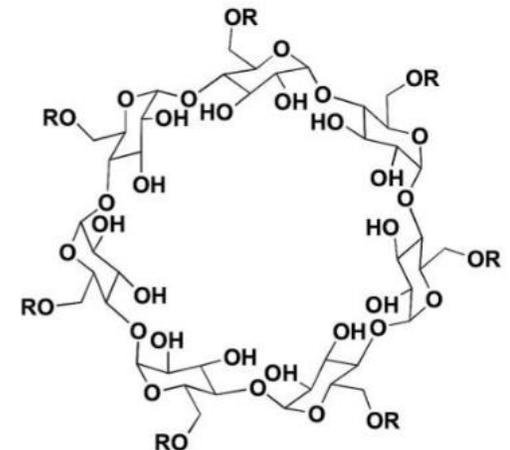
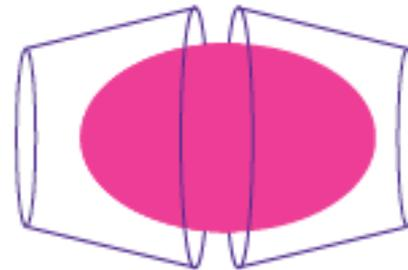


## Cyclodextrins strongly gained importance during past years

- Cyclodextrins are cyclic oligosaccharides derived from natural starch
- Units arranged in the form of hollow cone with hydrophilic exterior and hydrophobic cavity
- First discovered in the 19<sup>th</sup> century; nowadays increasing use as pharmaceutical excipient

## Benefits of Cyclodextrins<sup>[1]</sup>:

- Increases solubility and stability
- Enhances bioavailability
- Masks odors and tastes
- Reduces aggregation behavior of biologics



[1] Cyclodextrins, Kurkov & Loftsson, 2012



## Derivates offer enhanced solubility

- Currently, there are >120 products formulated with Cyclodextrins on the market
- SBE- $\beta$ -CD containing drugs nearly doubled during past years

Route	$\alpha$ -CD	$\beta$ -CD	$\gamma$ -CD	RM- $\beta$ -CD	HP- $\beta$ -CD	SBE- $\beta$ -CD	HP- $\gamma$ -CD
Oral	3 [0]	37 [0]	2 [0]	1 [0]	12 [1]	0 [1]	2 [0]
Buccal	0 [0]	0 [0]	0 [0]	0 [0]	4 [0]	0 [0]	0 [0]
Topical	1 [0]	7 [0]	0 [0]	0 [0]	1 [0]	0 [0]	0 [0]
Ophthalmic	0 [0]	1 [0]	0 [0]	1 [0]	2 [0]	0 [0]	0 [0]
Parenteral	6 [0]	0 [0]	0 [0]	0 [0]	19 [3]	25 [5]	1 [0]
<b>Total</b>	<b>10 [0]</b>	<b>45 [0]</b>	<b>2 [0]</b>	<b>2 [1]</b>	<b>38 [4]</b>	<b>25 [6]</b>	<b>3 [0]</b>

Marketed products employing the respective CD; numbers in brackets indicate new products currently in clinical phase  
 Not shown here: studies ongoing to broaden indication of existing drugs; PharmaCircle database, March 2022

RM: randomly methylated; HP: hydroxypropyl; SBE: sulfobutylether



## HP- $\beta$ -CD and SBE- $\beta$ -CD have very good safety profiles

- Part of FDA list of inactive ingredients
- Marketed drugs exist for parenteral administration
- Usually renally excreted intact without metabolization<sup>[1]</sup>
- Shielding effects can help minimizing gastrointestinal irritation<sup>[1]</sup>
- Very versatile for liquid, semi-solid and solid formulations

## Cyclodextrins and COVID-19

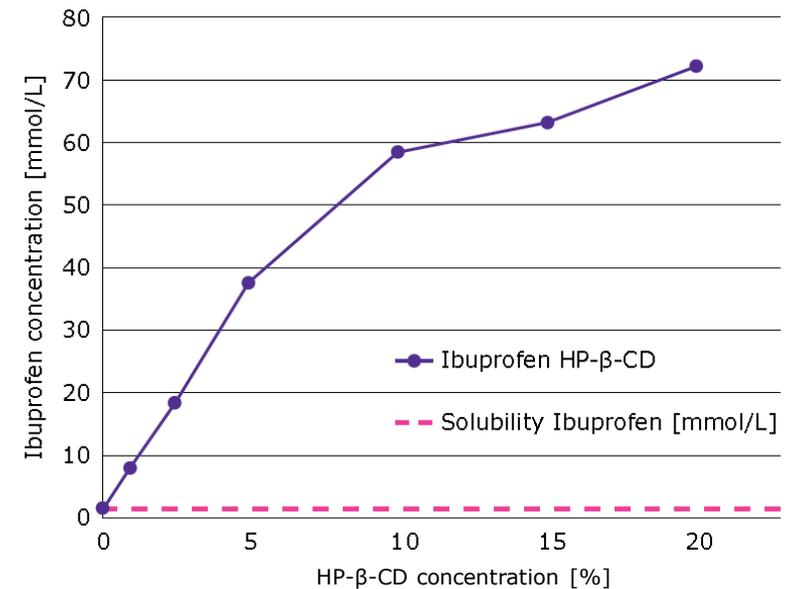
- Remdesivir (Gilead): SBE- $\beta$ -CD to increase solubility<sup>[2]</sup>
- COVID-19 vaccine (J&J): HP- $\beta$ -CD to protect proteins<sup>[3]</sup>

HP: hydroxypropyl; SBE: sulfobutylether

[1] Questions and answers on cyclodextrins used as excipients in medicinal products for human use, EMA, 2017

[2] Sulfobutylether- $\beta$ -cyclodextrin, Stella & Rajewski, 2020

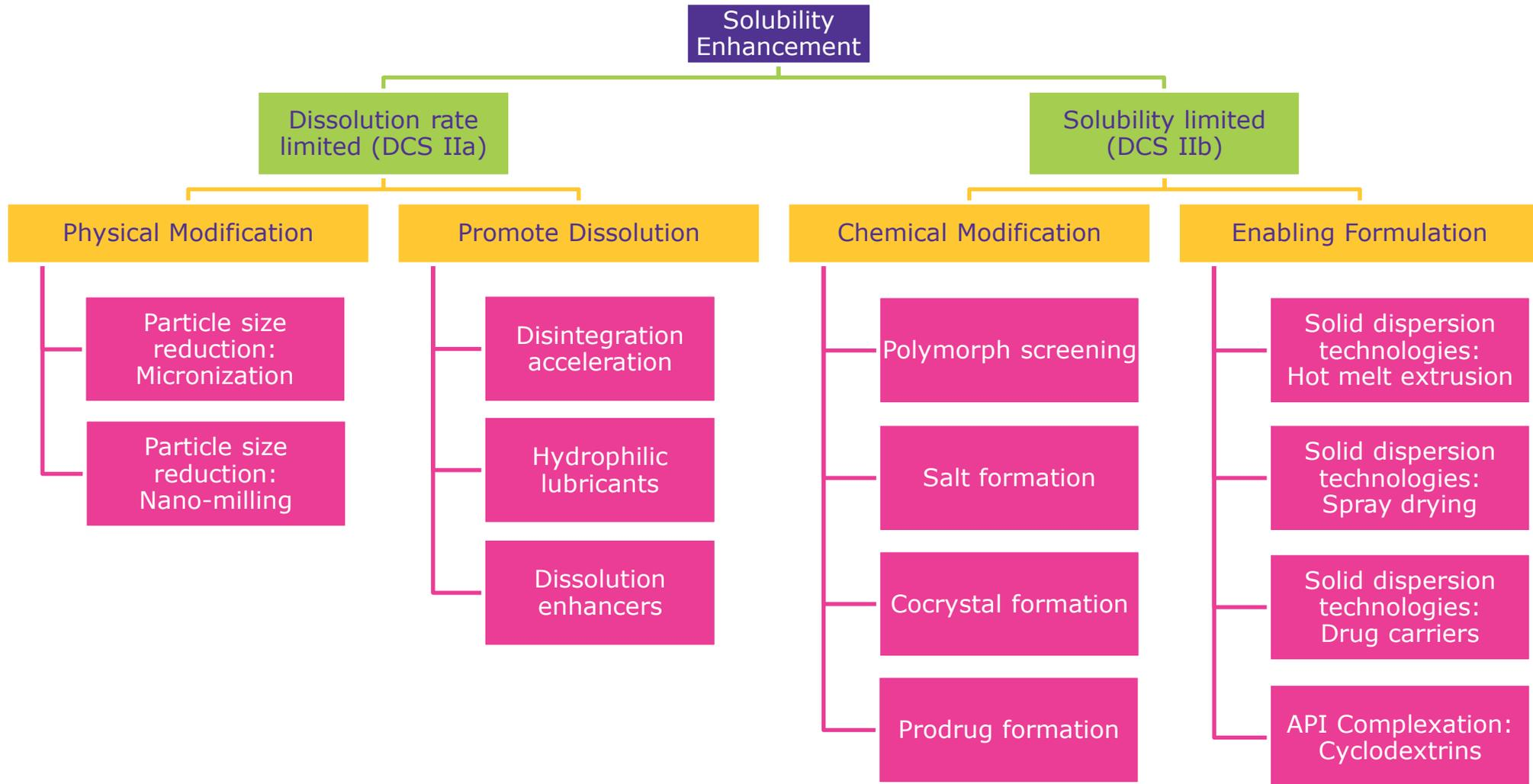
[3] Cyclodextrins in Antiviral Therapeutics and Vaccines, Braga *et al.*, 2021



Solubility of Ibuprofen as a function of HP- $\beta$ -CD concentration, internal data



# Strategies and Enabling Technologies for Enhancing API Solubility



**SAFC**®

Pharma & Biopharma  
Raw Material Solutions

**Dr. Thomas Briel**

Strategic Marketing Manager –  
Excipients Liquid Application

*thomas.briel@emdgroup.com*

**Dr. Markus Lubda**

Strategic Marketing Manager -  
Excipients Solid Application

*markus.lubda@emdgroup.com*

MilliporeSigma, the Vibrant M, Parateck and SAFC are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners.

