

**Sigma-Aldrich®**

Lab & Production Materials

**MERCK**



# Faster discovery Faster recovery

High-Throughput Covidytm™  
Substrates & ACE2 Kits for  
Covid-19 Drug Discovery

The right assay kit for your metabolite  
and Enzyme Analysis

We provide a comprehensive portfolio  
of assay kits for a broad range of  
applications in life science research,  
drug discovery and development, and  
environmental monitoring and testing.

- easy-to-use, precise, and highly reliable assays
- spectrophotometric and/or fluorometric detection methods
- also suitable for high-throughput analysis
- Each lot is tested for suitability, reproducibility and quality to ensure



As Covid-19 continues to impact the world, scientists are rushing to lives. We are committed to helping you reach your goals faster. Our extensive range of metabolomics products now includes high-quality Covidytm™ substrates and ACE2 kits that allow high-throughput handling for more rapid and reliable Discover our unparalleled portfolio for your breakthroughs in coronavirus therapy.

## Sample Preparation

Kit #	Product
MAK191	Carrez Clarification Reagent Kit
MAK227	Chloramphenicol Extraction Kit
MAK175	Cholesterol Extraction Kit
MAK341	Deproteinizing Sample Preparation Kit
MAK342	Deproteinizing Sample Preparation Kit II
MAK226	Enrofloxacin Extraction Kit
MAK372	Fatty Acid Extraction Kit for Pet Food
MAK338	Fatty Acid Extraction Kit, High Standard (450 ug/ml)

Kit #	Product
MAK174	Fatty Acid Extraction Kit
MAK225	Fatty Acid Methyl Ester Purification Kit Manufactured by Nacalai Tesque, Inc.
MAK224	Fatty Acid Methylation Kit Manufactured by Nacalai Tesque, Inc.
MAK339	Lipopolysaccharide (LPS) Isolation Kit
MAK340	Microsome Isolation Kit
MAK343	PEG Virus Precipitation Kit

## Covidytm™ Substrates

Covidytm™ are peptide substrates with a specific amino acid sequence that can be cleaved by coronavirus proteases. The dark-FRET peptide contains a quencher and a fluorescent donor on the N- and C-terminals, respectively, where the fluorescence of the donor is effectively quenched when the peptide is intact. However, when the peptide is hydrolyzed by coronavirus proteases, the fluorescent donor fragment generates significantly enhanced fluorescence. In this way, activity of coronavirus proteases can be effectively monitored through the fluorescence intensity of the fluorescent donor.

- Suitable for developing methods to screen potential inhibitors of coronavirus proteases
- Compatible with high-throughput handling systems

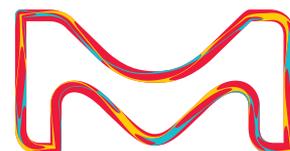
Order No.	Product Name	Description
MAK373	Covidytm™ EN450 Substrate	Peptide substrate containing a 14 amino acid sequence that can be cleaved by coronavirus proteases
MAK374	Covidytm™ ED450 Substrate	Peptide substrate containing a 12 amino acid sequence that can be cleaved by coronavirus proteases
MAK375	Covidytm™ TF670 Substrate	Peptide substrate containing a 14 amino acid sequence that can be cleaved by coronavirus proteases
MAK376	Covidytm™ IF670 Substrate	Peptide substrate containing a 12 amino acid sequence that can be cleaved by coronavirus proteases

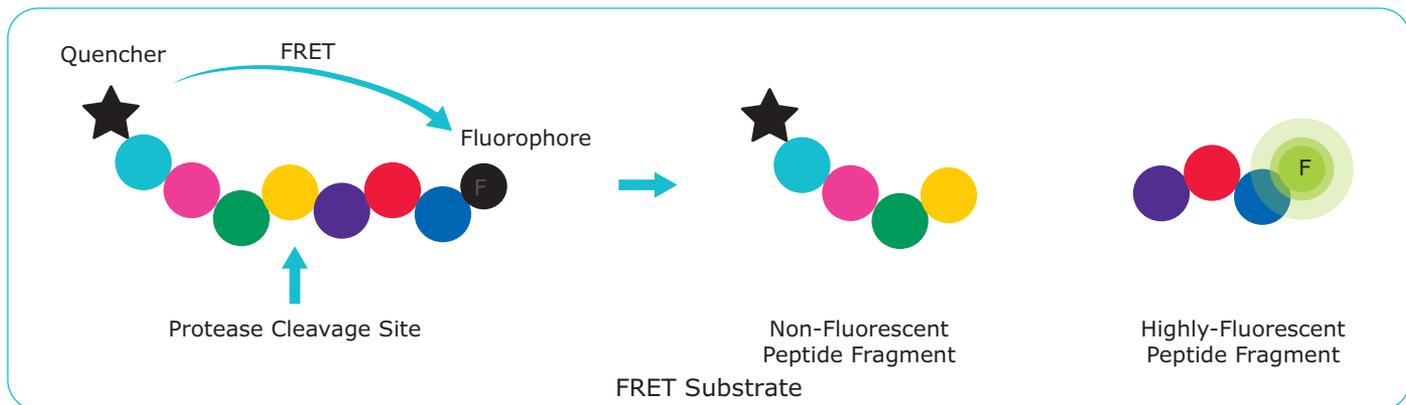
## ACE2 Activity Assay Kit & Inhibitor Screening Kit

The Angiotensin II Converting Enzyme (ACE2) Activity Assay Kit and Screening Kit both use the ability of an active ACE2 to cleave a synthetic MCA-based peptide substrate to release a free fluorophore. Released MCA can be easily quantified using a fluorescence microplate reader. The kits also include an ACE2-specific inhibitor that can differentiate ACE2 activity from other proteolytic activity.

- Activity Assay Kit: suitable for detection of ACE2 activity in tissue (e.g. lung, heart and kidney) and overexpressed ACE2 in cell lysates, as well as for determination of enzymatic activity of purified ACE2
- Inhibitor Screening Kit: suitable for screening/characterizing inhibitors/ligands of ACE2
- Both kits are compatible with high-throughput handling systems

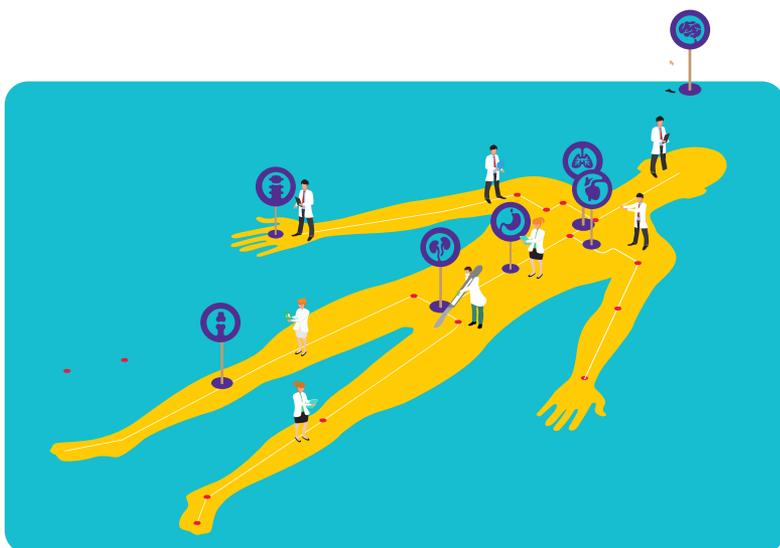
Order No.	Product Name	Description
MAK377	Angiotensin II Converting Enzyme (ACE2) Activity Assay Kit	Easy-to-use kit for detection of ACE2 concentrations as low as 0.4 mU
MAK378	Angiotensin II Converting Enzyme (ACE2) Inhibitor Screening Kit	Convenient kit used to screen for potent inhibitors of ACE2 activity





More products available at: [SigmaAldrich.com/covid-19](https://www.sigmaaldrich.com/covid-19)

To learn more about additional metabolism assay kits for your drug discovery research, visit [SigmaAldrich.com/metabolism-assay-kits](https://www.sigmaaldrich.com/metabolism-assay-kits)



### Contact Us

PT Merck Chemicals and Life Sciences  
 Arkadia Green Park - G building, 9th floor  
 Jl. TB Simatupang No.Kav. 88, Pasar Minggu Jakarta, 12520, Indonesia  
[www.merckmillipore.com](http://www.merckmillipore.com) | [www.sigmaaldrich.com](http://www.sigmaaldrich.com)

PIC : Sri Hayuni (HP:0811-8066-201)

Email : [sri.hayuni@merckgroup.com](mailto:sri.hayuni@merckgroup.com)

Toll Free: 0800 140 1253 | [CS.Merck@merckgroup.com](mailto:CS.Merck@merckgroup.com)

[@mercklifescience.id](https://www.instagram.com/mercklifescience.id) | [Merck Life Science Indonesia](https://www.youtube.com/channel/UC...)

[@merckindonesia](https://www.line.me/@merckindonesia) | [Merck Life Science Indonesia](https://www.linkedin.com/company/merck-life-science-indonesia)

© 2021 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck, the vibrant M, and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

