

Product Information

KitAlysis™ High-Throughput Metal Scavenger Reaction Screening Kit

Catalog Number **KITALYSIS-SCAV**

Storage Temperature 20–25 °C

TECHNICAL BULLETIN

Product Description

The KitAlysis™ High-Throughput Metal Scavenger Reaction Screening Kit was designed for quick selection of scavenger in removing excess metals from reactions or isolated product. Specifically chosen by chemistry experts, each of the twelve scavengers is provided in 10 mg and 50 mg amounts to quickly assess not only the optimal scavenger, but the optimal w/w% of scavenger to substrate as well.

Each Metal Scavenger Reaction Screening Kit contains 2 individual screening sets. Each screening set has components to run 24 unique metal scavenger reactions useful for:

- Broad-spectrum metal removal
- Impurity removal
- Color treatment

Access www.sigmaaldrich.com/kitalysis for a detailed step-by-step protocol.

Screening Kit Components

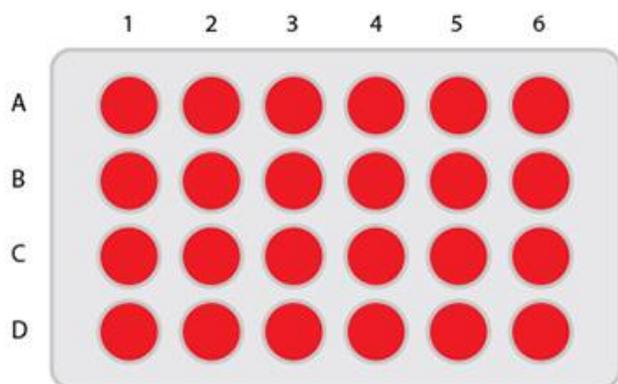
Two screening sets individually sealed under nitrogen in mylar bags. Each bag includes:

- 12 × 2 pre-weighed scavengers/resins in glass vials topped with cap mat, see Figures 1 and 2.

Figure 1.
Scavengers

Name	Catalog Number	Vials
Cysteine functionalized Silica Gel	900677	A1, C1
<i>N</i> -Propyldiethanolamine-functionalized silica gel	900676	A2, C2
Triaminetetraacetate, Na salt-functionalized silica gel	900674	A3, C3
Triaminetetraacetic acid-functionalized silica gel	900675	A4, C4
DMT-functionalized silica gel	900680	A5, C5
3-(Diethylenetriamino)propyl-functionalized silica gel	537926	A6, C6
Mixture of metal scavengers on silica gel	900673	B1, D1
DOTA-functionalized silica gel	900679	B2, D2
AMPA-functionalized silica gel	900678	B3, D3
3-(Ethylenediamino)propyl-functionalized silica gel	537918	B4, D4
3-(1-Thioureido)propyl-functionalized silica gel	569879	B5, D5
3-Mercaptopropyl-functionalized silica gel	538086	B6, C6

Figure 2.
Position of Scavengers



Reagents and Equipment Required, but Not Provided.

- 1 mL syringe with needle or 1 mL pipette with tips
- 50 mL flask
- Solvent [DCM, THF, EtOAc, DMF, MeOH, Toluene, Water (pH 2-10) or any other solvent in which the substrate is soluble]
- Substrate: metal contaminated crude or purified material
- Orbital shaker for 96 well plates (the use of stir bar is not recommended with some of the resins)

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

Store the High-Throughput Metal Scavenger Reaction Screening Kit at room temperature.

KitAlysis is a trademark of Sigma-Aldrich Co. LLC.

DR,MAM 10/17-1