

Succinic acid disodium salt

Succinic acid is a dicarboxylic acid that plays a role in the citric acid cycle, yielding energy. It is used as an acidity regulator in the food and beverage industry, and sold as a food additive and supplement.

Materials Provided

Catalog Number:	92-1193
Quantity Supplied:	1 vial(s), 10 mg (re-packaged) per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	162.05 g/mol
Source:	Synthetic
Purity:	≥ 98%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	RT Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	25 mM
Reconstitution Volume:	2.468 mL
Reconstitution Solvent:	H ₂ O

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com