

α -Ketoglutaric acid sodium salt

α -Ketoglutaric acid is a naturally occurring ketone derivative of glutaric acid and intermediate of the krebs cycle. It serves as a nitrogen transporter, antioxidant, anabolic, wound healer, and may have anti-aging properties. Used to treat kidney disease, GI disorders, bacterial infections, liver problems, cataracts, and yeast infections. Commonly sold as a supplement to help combat ammonia toxicity.

Materials Provided

Catalog Number:	92-1212
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:	168.08 g/mol
Source:	Synthetic
Purity:	$\geq 98\%$
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	2-8°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	100 mM
Reconstitution Volume:	0.595 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