

3-Hydroxyoctanoic acid

3-Hydroxyoctanoic acid is a signaling chemical emitted by orchids that may also be a marker for primary defects of the β -hydroxy fatty acid metabolism. A polymer of 3-hydroxyoctanoic acid produces a carbon and energy reserve form for bacteria. Also used to produce polyhydroxyalkanoates used in medical devices and tissue engineering applications.

Materials Provided

Catalog Number:	92-1189
Quantity Supplied:	1 vial(s), 5 mg 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:	160.21 g/mol
Source:	Synthetic
Purity:	$\geq 96.5\%$ by GC
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:	50 mM
Reconstitution Volume:	0.624 mL
Reconstitution Solvent:	Ethanol

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