

Triiodothyronine (T3)

Triiodothyronine (T3) is a thyroid hormone that is an active form of thyroxine. It affects nearly every physiological process in the body, and has a vital role in metabolism, heart and digestive function, muscle control, brain development, and bone maintenance. Agonist of the thyroid hormone receptors.

Materials Provided

Catalog Number:	92-1137
Quantity Supplied:	1 vial(s), 10 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:	672.96 g/mol
Source:	Synthetic
Purity:	≥ 95%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	5 mM
Reconstitution Volume:	2.972 mL
Reconstitution Solvent:	1N NaOH

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