

## Uridine 5'-Triphosphate

Uridine 5'-Triphosphate (UTP) is a uracil nucleoside triphosphate that acts as a substrate for RNA synthesis during transcription. UTP also provides a source of energy and activates substrates in metabolic reactions, like ATP.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1109</b>
<b>Quantity Supplied:</b>	1 vial(s), 10 mg per vial
<b>Reconstitution Buffer:</b>	1 mL

### Description

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

### Product Information

<b>Molecular Weight:</b>	484.14 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 90%
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	≤ -20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	50 mM
<b>Reconstitution Volume:</b>	0.413 mL
<b>Reconstitution Solvent:</b>	Reconstitution Buffer

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)