

## Serotonin 5-HT

Serotonin, or 5-hydroxytryptamine (5-HT), is a mono-amine neurotransmitter derived from tryptophan that is responsible for regulation of mood, commonly referred to as the "depression gene". It is primarily located in the CNS and GI tract, and modulates physiological functions related to neurotransmission, GI motility, pain perception, hemostasis, hunger, and cardiovascular integrity.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1108</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>	212.68 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 98%
<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>	0.94 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)