

## LCK Inhibitor

LCK (lymphocyte-specific protein tyrosine kinase) inhibitor is a pyrrolopyrimidine that block the activity of two forms of Lck kinase, Lck (64-509) and Lckcd, and has been shown to inhibit T cell receptor-stimulated IL-2 production. The LCK protein is a member of the Src-family of non-receptor protein tyrosine kinases and plays a critical role in the initial steps of T cell receptor signaling that trigger the production of cytokines.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1322</b>
<b>Quantity Supplied:</b>	1 vial(s), 1 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

<b>Molecular Weight:</b>	370.5
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥95% by HPLC
<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>	10 mM
<b>Reconstitution Volume:</b>	0.270 mL
<b>Reconstitution Solvent:</b>	DMSO

***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)