

Recombinant Human IL-2, Cell Culture Grade

Interleukin 2 (IL-2) is a cytokine protein and intricate member of the immune response. Regulates activation of T and B lymphocyte proliferation, and also stimulates monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells. This cell culture grade recombinant human IL-2 is a variant form of IL-2 (contains a serine to cysteine substitution) that can be used for differentiation studies, functional assays, and cell culture.

Materials Provided

Catalog Number:	92-1331
Quantity Supplied:	1 vial(s), 10 µg per vial
Reconstitution Buffer:	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

Molecular Weight:	15,400 g/mol
Source:	Recombinant
Purity:	>97%
Endotoxin Level:	<0.10 EU
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:	100 µg/mL
Reconstitution Volume:	0.100 ml
Reconstitution Solvent:	Reconstitution Buffer

Handling Instructions:

specific activity of $\geq 5.0 \times 10^6$ IU/mg

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