

## Recombinant Human SCF

Stem cell factor (SCF) is a cytokine with important roles in hematopoiesis, spermatogenesis, and melanogenesis. It potently binds the c-KIT receptor (CD117) and other cytokines to preserve the viability and facilitate the proliferation and differentiation of hematopoietic stem and progenitor cells.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1114</b>
<b>Quantity Supplied:</b>	1 vial(s), 2 µg 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>	18400 g/mol
<b>Source:</b>	Recombinant
<b>Purity:</b>	≥ 98% by HPLC
<b>Endotoxin Level:</b>	< 0.1 ng/µg (<1EU/µg).
<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 µg/mL
<b>Reconstitution Volume:</b>	0.040 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)