

Stem cell factor (human), (recombinant)

Stem cell factor (SCF) is a cytokine made by fibroblasts and endothelial cells. SCF binds to the receptor known as c-Kit (CD117) and is thought to play a critical role in the maintenance or survival of hematopoietic stem cells. Human SCF shows no activity on mouse cells, but mouse and rat SCF are active on human cells.

Ordering Information

[Order Online »](#)

ALX-201-804-0010	10µg
ALX-201-804-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

- High purity
- Carrier-free

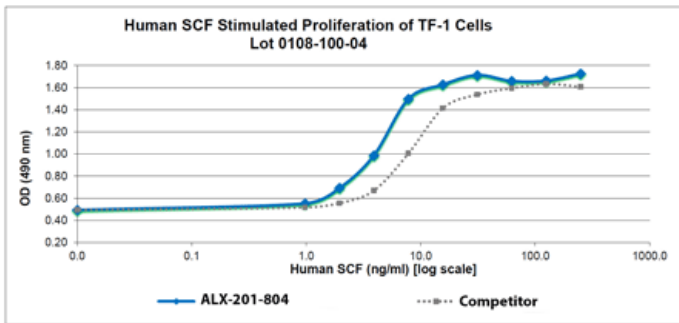


Figure 1: Human SCF Bioactivity Data. Serial dilutions of human SCF, starting at 250ng/ml, were added to TF-1 cells. After 68 hours, cell proliferation was measured and the linear portion of the curve was used to calculate the ED50. The ED50 for the lot shown of human SCF is 3.3-5ng/ml. This value is comparable to the typical expected value of 1-5ng/ml.

Handling & Storage

Use/Stability	It is recommended that a carrier protein (0.1% HSA or BSA) is added for long term storage.
Handling	Centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	SCF, c-Kit ligand, KL, Steel factor, MGF
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the dose-dependent stimulation of human TF-1 cells which is typically observed at a concentration of ≤ 15 ng/mL.
Endotoxin Content	≤ 1 EU/ μ g protein measured by kinetic LAL analysis.
Formulation	Lyophilized from a 0.2 μ m filtered solution containing 10 mM Sodium Phosphate, 50 mM Sodium Chloride, pH 7.5.
MW	~18.4kDa
Purity	$\geq 97\%$ (SDS-PAGE using Coomassie Staining)
Reconstitution	When reconstituting the product, gently pipet and wash down the sides of the vial to ensure full recovery of the protein into solution. It is recommended to reconstitute the lyophilized product with sterile water at a concentration of 0.1 mg/ml, which can be further diluted into other aqueous solutions.
Source	Produced in <i>E. coli</i> . Non-glycosylated protein, containing 165 amino acids.
UniProt ID	P21583

Last modified: August 6, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)