

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-201-804-0002 | 2µg |
| ALX-201-804-0010 | 10µg |
| ALX-201-804-0100 | 100µg |

Manuals, SDS & CofA

[View Online »](#)

- High purity
- Carrier-free

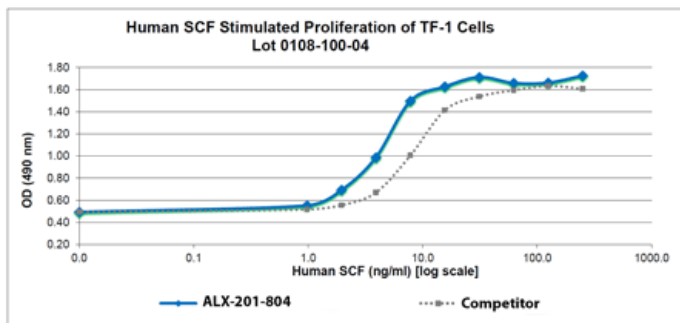


Figure 1: Human SCF Bioactivity Data. Serial dilutions of human SCF, starting at 250ng/ml, were added to TF-1 cells. After 68 hours, cell proliferation was measured and the linear portion of the curve was used to calculate the ED50. The ED50 for the lot shown of human SCF is 3.3-5ng/ml. This value is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



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)