

Stem cell factor (mouse), (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-201-805-0002 | 2µg |
| ALX-201-805-0010 | 10µg |
| ALX-201-805-0100 | 100µg |

Manuals, SDS & CofA

[View Online »](#)

- High purity
- Carrier-free

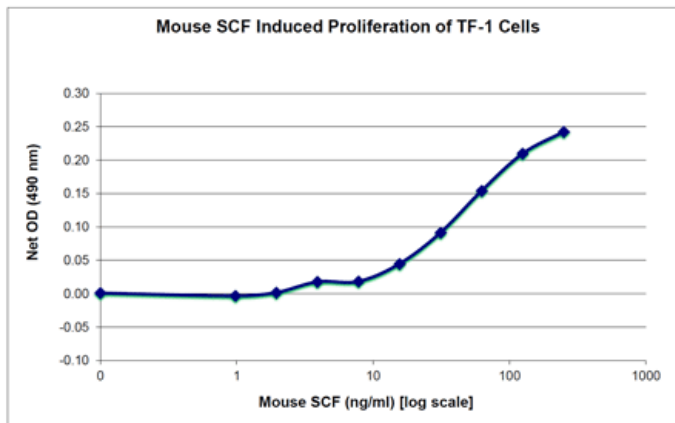


Figure 1: Mouse SCF Bioactivity Data. Serial dilutions of mouse SCF, starting at 250ng/ml, were added to TF-1 cells growing in GM-SCF free media. Cell proliferation was measure after 65 hours and the linear portion of the curve was us used to calculate the ED50. The ED50 for this lot of mouse SCF was between 34-51ng/ml. The typical range is between 2.5-10ng/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 proliferation of human TF-1 cells which is typically observed at concentrations <10ng/ml. |
| Endotoxin Content | ≤1 EU/μg protein measured by kinetic LAL analysis. |
| Formulation | Lyophilized from a sterile aqueous solution containing 20mM sodium phosphate, pH 6.5. |
| MW | ~18.3kDa |
| Purity | ≥95% (reducing and non-reducing SDS-PAGE) |
| 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 164 amino acids. |
| UniProt ID | P20826 |



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)