

Flt-3 ligand (human), (recombinant)

FMS-related tyrosine kinase 3 ligand (Flt-3 ligand) is a growth factor important for the proliferation of hematopoietic cells. Flt-3 ligand binds to, and transmits signals through, the receptor tyrosine kinase known as FMS-like tyrosine kinase-3 (Flt-3). Flt-3 ligand promotes long-term expansion and differentiation of human pro-B cells in the presence of IL-7 or in combination of IL-7 and IL-3. Human Flt-3 ligand can stimulate the proliferation of cells expressing mouse Flt-3 receptors.

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

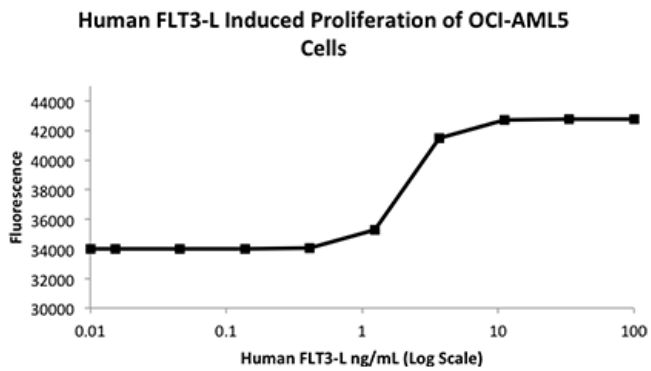


Figure 2: Serial dilutions of Human FLT-3 Ligand, starting at 100 ng/mL, were added to OCI-AML5 cells. After 60 hours, proliferation can be measured and the linear portion of the curve can be used to calculate the ED50.

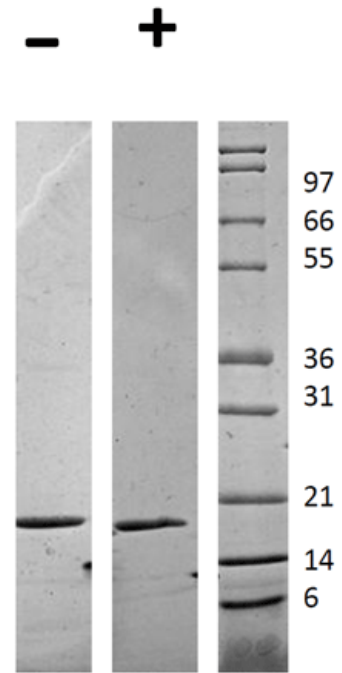


Figure 1: 1 μ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human Flt-3 ligand has a predicted MW of 17.6 kDa.

Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | It is recommended that a carrier protein (0.1% 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 | FMS-related tyrosine kinase 3 ligand |
| Appearance | White lyophilized (freeze-dried) powder. |
| Biological Activity | The activity is determined by the dose-dependent stimulation of the proliferation of human Aml5 cells which is typically observed at a concentration of 1-3 ng/mL. |
| Endotoxin Content | ≤1 EU/μg protein measured by kinetic LAL analysis. |
| Formulation | Lyophilized from sterile filtered solution containing 10 mM sodium phosphate, 50 mM NaCl, pH 7.5. |
| MW | ~17.6kDa |
| Purity | ≥95% (Reducing and Non-reducing SDS-PAGE) |
| Reconstitution | Reconstitute with sterile water at 0.1 mg/ml. |
| Source | Produced in <i>E. coli</i> . Non-glycosylated protein, containing 155 amino acids. |
| UniProt ID | P49771 |