

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.

Ordering Information

[Order Online »](#)

ALX-201-800-0002	2µg
ALX-201-800-0010	10µg
ALX-201-800-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

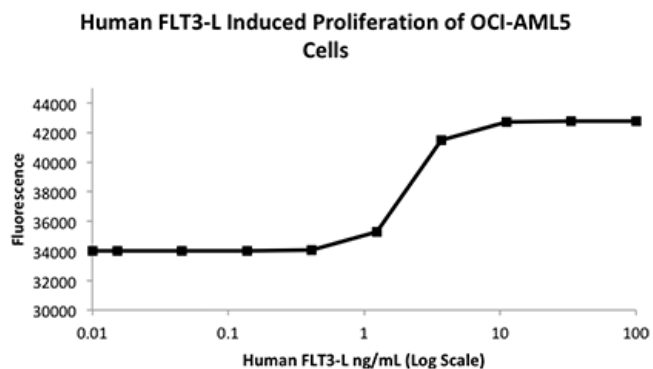


Figure 2: Serial dilutions of Human FLT-3 Ligand, starting at 100 ng/mL, were added to OCI-AML5 cells. After 60 hours, proliferation can be measured and the linear portion of the curve can be used to calculate the ED50.

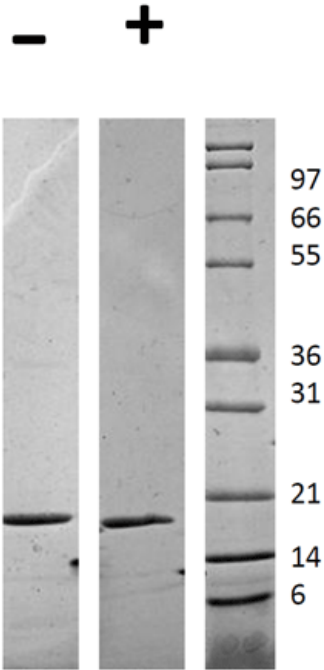


Figure 1: 1µg in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human Flt-3 ligand has a predicted MW of 17.6 kDa.

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