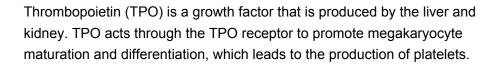
Thrombopoietin (mouse), (recombinant)



• Carrier-free

Ordering Information

Order Online »

ALX-201-808-0002	2μg
ALX-201-808-0010	10µg

Manuals, SDS & CofA

View Online »

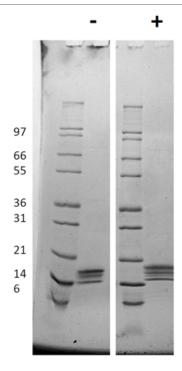


Figure 2: 1 μ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Mouse TPO has a predicted MW of 18.7 kDa.

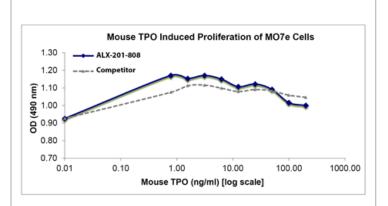


Figure 1: Mouse TPO Bioactivity Data. Serial dilutions of mouse TPO, starting at 200ng/ml, were added to MO7e cells growing in the presence of 1ng/ml IL-3 and 2.5ng/ml SCF. Proliferation was measure after 5 days and the linear portion of the curve was us used to calculate the ED50. The ED50 for this lot of mouse TPO was less than 0.8ng/ml. This value is comparable to the typical expected range of 1ng/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 TPO, Megakaryocyte colony stimulating factor, c-MPL

ligand, MGDF

Appearance White lyophilized (freeze-dried) powder.

Biological Activity The activity is determined by the dose-dependent

stimulation of MO7e cells which is typically observed at

concentations <1ng/ml.

Endotoxin Content ≤1 EU/µg protein measured by kinetic LAL analysis.

Formulation Lyophilized. Sterile filtered.

MW ~18.7kDa

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 *E. coli*. Non-glycosylated protein, containing

174 amino acids (which comprise the receptor binding

domain).

UniProt ID P40226

Last modified: May 29, 2024

