

Thrombopoietin (mouse), (recombinant)

Thrombopoietin (TPO) is a growth factor that is produced by the liver and kidney. TPO acts through the TPO receptor to promote megakaryocyte maturation and differentiation, which leads to the production of platelets.

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

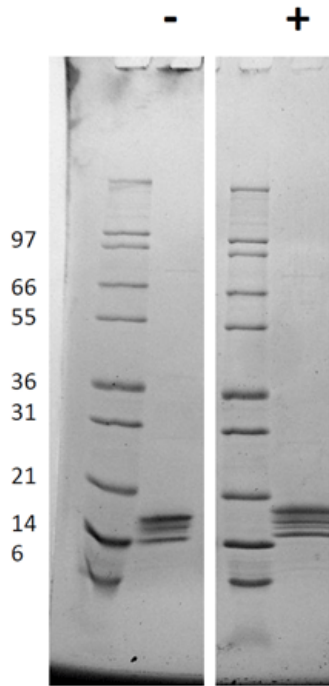


Figure 2: 1 μ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Mouse TPO has a predicted MW of 18.7 kDa.

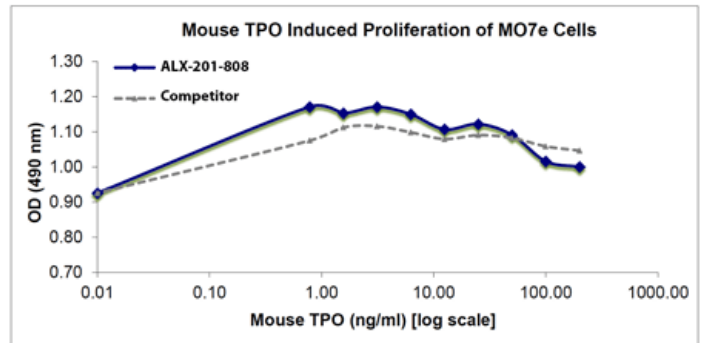


Figure 1: Mouse TPO Bioactivity Data. Serial dilutions of mouse TPO, starting at 200ng/ml, were added to MO7e cells growing in the presence of 1ng/ml IL-3 and 2.5ng/ml SCF. Proliferation was measure after 5 days and the linear portion of the curve was us used to calculate the ED50. The ED50 for this 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 concentrations <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 <i>E. coli</i> . Non-glycosylated protein, containing 174 amino acids (which comprise the receptor binding domain).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com