

# LIF (mouse), (recombinant)

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-201-242-L001	1ml
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** Stable for 6 months when stored at +4°C.

**Handling** Avoid freeze/thaw cycles.

**Short Term Storage** +4°C

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Leukemia inhibitory factor

**Biological Activity** Induces macrophage differentiation and proliferation of hematopoietic stem cells; maintains undifferentiated state of embryonic stem (ES) cell cultures; inhibits proliferation of the mouse myeloid cell line M1.

**Formulation** Liquid. 0.2µm-filtered solution in PBS containing 0.5% BSA. Contains no preservatives. Further dilutions should be made with medium containing 1% BSA.

**MW** ~20kDa

**Purity** ≥95% (SDS-PAGE)

**Purity Detail** Purified by FPLC.

**Quantity** 10<sup>7</sup>U

**Source** Produced in *E. coli*.

**Technical Info / Product Notes** **Maintenance of ES Cells:** 1,000U/ml are recommended to maintain undifferentiated ES cells, R1 or E14 cells.

**M1Growth Inhibition Assay:** 50units/ml are defined as the concentration of LIF that induces 50% of differentiated cells in tissue culture medium.



ENZO LIFE SCIENCES  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

P09056

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)