

# 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 <i>E. coli</i> .

Technical Info / Product Notes	<p><b>Maintenance of ES Cells:</b> 1,000U/ml are recommended to maintain undifferentiated ES cells, R1 or E14 cells.</p> <p><b>M1Growth Inhibition Assay:</b>50units/ml are defined as the concentration of LIF that induces 50% of differentiated cells in tissue culture medium.</p>
--------------------------------	--

Last modified: May 29, 2024



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

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)