

# Fc (human):Dectin-2 (mouse), (recombinant)

Dectin-2 (dendritic cell-associated C-type lectin-2) is a lectin that has been reported to bind to zymosan as well as to hyphae of organisms such as *Candida albicans*. Dectin-2 exhibits specific recognition of high-mannose structures. Activation of Dectin-2 leads to phosphorylation of FcRg and up-regulated secretion of TNF- $\alpha$  and IL-1ra. Dectin-2 might act on inflammatory macrophages and DC to detect pathogenic organisms and possibly endogenous glycoproteins.

## Ordering Information

[Order Online »](#)

ALX-522-126-C050	50 $\mu$ g
------------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 6 months after receipt when stored at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Dendritic cell-associated lectin 2
<b>Concentration</b>	1mg/ml after reconstitution.
<b>Endotoxin Content</b>	<0.1EU/μg purified protein (LAL test; Associates of CAPE COD Inc.).
<b>Formulation</b>	Lyophilized. Contains PBS.
<b>MW</b>	~55kDa (SDS-PAGE)
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute with 50μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.
<b>Source</b>	Produced in HEK 293 cells. The extracellular domain of dectin-2 (aa 44-209) is fused at the N-terminus to the Fc portion of human IgG1.
<b>UniProt ID</b>	Q9JKF4



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-  
fr@enzolifesciences.com](mailto:info-fr@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)