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 upregulated secretion of TNF- α 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

50µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 6 months after receipt when stored at -20°C.

Handling Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dendritic cell-associated lectin 2

Concentration 1mg/ml after reconstitution.

Endotoxin Content <0.1EU/µg purified protein (LAL test; Associates of CAPE

COD Inc.).

Formulation Lyophilized. Contains PBS.

MW ~55kDa (SDS-PAGE)

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute with 50µl sterile water. Further dilutions

should be made with medium containing 5% fetal calf

serum.

Source 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.

UniProt ID Q9JKF4

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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