IL-35 (human):Fc (human), (recombinant)

Highly pure interleukin involved in T cell regulation.

EBI3 (Epstein-Barr virus-induced gene 3 protein; EBV-induced gene 3 protein; IL-27B; IL-27 subunit β) and IL-12A (IL-12 subunit p35; CLMF p35) form the heterodimeric cytokine IL-35 (interleukin-35) involved in T cell regulation. IL-35 is highly expressed in regulatory T cell (Treg) but not in effector or naive T cells. IL-35 is required for the suppressive function of T-reg *in vitro* and *in vivo*.

Citations: 11

View Online »

Ordering Information

Order Online »

ALX-522-140-C010

10µg

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.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-35, EBI3:IL-12A, IL-27B:CLMF p35

Concentration 0.1mg/ml after reconstitution.

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

Cod).

Formulation Lyophilized. Contains PBS.

MW ~85kDa

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute with 100µl sterile water. Further dilutions

should be made with medium containing 5% fetal calf

serum or a carrier protein.

Source Produced in HEK 293 cells. Human EBI3 (aa 21-229)

linked to human IL-12A (aa 23-219) via the linker

GGGSGGGSGGGSGGS is fused at the C-terminus to

the Fc portion of human IgG1 and a FLAG®-tag.

Technical Info / Product NotesHistorical data has shown that IL-35 binds to IL-12Rß2,

and it has been demonstrated in the literature to inhibit

mononuclear cell proliferation.

FLAG is a registered trademark of Sigma-Aldrich Co.

Last modified: May 29, 2024



info-

ENZO LIFE SCIENCES, European Sales Office ENZO LIFE SCIENCES Phone: 800.942.0430 (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com France Phone: +33 472 440 655 fr@enzolifesciences.com

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

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