

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	10 μ g
------------------	------------

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 GGGSGGGSGGGSGGGS is fused at the C-terminus to the Fc portion of human IgG1 and a FLAG [®] -tag.
Technical Info / Product Notes	Historical data has shown that IL-35 binds to IL-12Rβ2, and it has been demonstrated in the literature to inhibit mononuclear cell proliferation. FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

Q14213 (IL-27B), P29459 (IL-21A)
European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

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

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