

PRODUCT INFORMATION



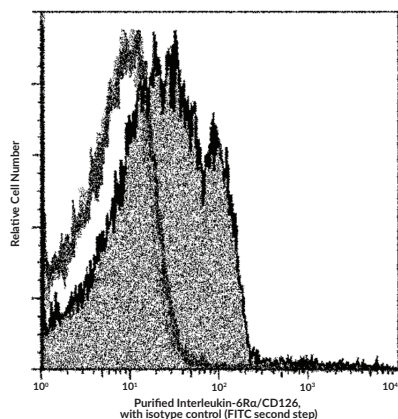
Interleukin-6R α /CD126 Rabbit Monoclonal Antibody (Clone 033)

Item No. 38089

Overview and Properties

Contents:	This vial contains 50, 100 μ l, or 1 ml of protein A-affinity purified monoclonal antibody
Synonyms:	IL-6R α , Interleukin-6 Receptor Subunit α
Immunogen:	Recombinant mouse IL-6R α
Cross Reactivity:	(+) IL-6R α , sIL-6R α
Species Reactivity:	(+) Mouse
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	0.2 μ m filtered solution in PBS
Clone:	033
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA (capture) and Flow cytometry (FC); the recommended starting dilution is 1:250-1:2,000 for ELISA (capture) and 1:25-1:100 for FC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Flow cytometric analysis of mouse Interleukin-6R α /CD126 expression on BABL/c splenocytes. Cells were stained with purified Interleukin-6R α /CD126 Rabbit Monoclonal Antibody (Clone 033), followed by a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/10/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Interleukin-6R α (IL-6R α), also known as CD126, is a type I cytokine receptor.¹ It is composed of an immunoglobulin-like domain, a cytokine-binding domain, a transmembrane domain, and a cytoplasmic domain and forms a heteromeric complex with IL-6 and gp130, which is the signal transducing subunit of the IL-6R that primarily acts through the JAK/STAT3 pathway.² IL-6R α is expressed in hepatocytes and monocytes and is localized to the plasma membrane but also exists as a soluble form (sIL-6R α) that is primarily generated *via* proteolytic cleavage of membrane-bound IL-6R α , but can also be formed *via* alternative splicing.³⁻⁵ Binding of IL-6 to membrane-bound IL-6R α , known as classical signaling, upregulates the synthesis of acute phase proteins, induces lipolysis in the liver, and induces the differentiation of T and B cells and is primarily anti-inflammatory.^{6,7} Trans-signaling consists of IL-6 binding to sIL-6R, which has various pro-inflammatory effects. Knockdown of *IL6R* increases sensitivity to tamoxifen in breast cancer cells.⁸ Formulations containing chimeric anti-IL-6R α monoclonal antibodies have been used in the treatment of rheumatoid arthritis, idiopathic arthritis, giant cell arteritis, cytokine release syndrome, and COVID-19. Cayman's Interleukin-6R α /CD126 Rabbit Monoclonal Antibody (Clone 033) can be used for ELISA (capture) and flow cytometry (FC) applications.

References

1. Hirano, T. Interleukin 6 and its receptor: Ten years later. *Int. Rev. Immunol.* **16(3-4)**, 249-284 (1998).
2. Salvati, A.L., Lahm, A., Paonessa, G., *et al.* Interleukin-6 (IL-6) antagonism by soluble IL-6 receptor α mutated in the predicted gp130-binding interface. *J. Biol. Chem.* **270(20)**, 12242-12249 (1995).
3. Bauer, J., Lengyel, G., Bauer, T.M., *et al.* Regulation of interleukin-6 receptor expression in human monocytes and hepatocytes. *FEBS Lett.* **249(1)**, 27-30 (1989).
4. Flynn, C.M., Kespohl, B., Daunke, T., *et al.* Interleukin-6 controls recycling and degradation, but not internalization of its receptors. *J. Biol. Chem.* **296**, 100434 (2021).
5. Riethmueller, S., Somasundaram, P., Ehlers, J.C., *et al.* Proteolytic origin of the soluble human IL-6R *in vivo* and a decisive role of N-glycosylation. *PLoS Biol.* **15(1)**, e2000080 (2017).
6. Del Giudice, M. and Gangestad, S.W. Rethinking IL-6 and CRP: Why they are more than inflammatory biomarkers, and why it matters. *Brain Behav. Immun.* **70**, 61-75 (2018).
7. Rose-John, S. Interleukin-6 family cytokines. *Cold Spring Harb. Perspect. Biol.* **10(2)**, a028415 (2018).
8. Tsoi, H., Man, E.P.S., Chau, K.M., *et al.* Targeting the IL-6/STAT3 signalling cascade to reverse tamoxifen resistance in estrogen receptor positive breast cancer. *Cancers (Basel)* **13(7)**, 1511 (2021).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM