

# PRODUCT INFORMATION

## Citrullinated Interleukin-6 (human, recombinant)

Item No. 30173

### Overview and Properties

**Synonyms:** B Cell Stimulatory Factor 2, BSF-2, CDF, CTL Differentiation Factor, HGF, Hybridoma Growth Factor, IFN- $\beta$ 2, IL-6, Interferon- $\beta$ 2

**Source:** Recombinant human N-terminal His-tagged IL-6 expressed in HEK293 cells, citrullinated by PAD4

**Amino Acids:** 30-212

**Uniprot No.:** P05231

**Molecular Weight:** 20.9 kDa

**Storage:** -80°C (as supplied)

**Stability:**  $\geq 1$  year

**Purity:** **batch specific** ( $\geq 85\%$  estimated by SDS-PAGE)

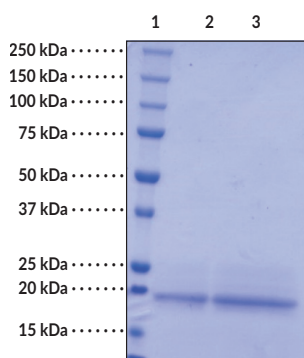
**Supplied in:** 50 mM HEPES, pH 8.0, with 150 mM sodium chloride, 10% glycerol, 2 mM  $\beta$ -mercaptoethanol

**Endotoxin Testing:**  $< 1.0$  EU/ $\mu$ g, determined by the LAL endotoxin assay

**Protein Concentration:** **batch specific** mg/ml

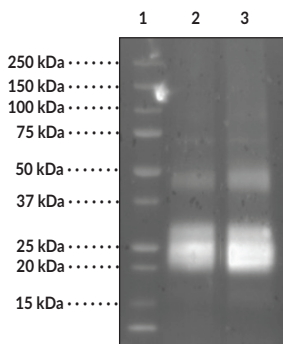
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: Citrullinated IL-6 (2  $\mu$ g)  
Lane 3: Citrullinated IL-6 (4  $\mu$ g)

SDS-PAGE Analysis of Citrullinated IL-6.



Lane 1: MW Markers  
Lane 2: Interleukin-6  
Lane 3: Citrullinated IL-6

Analysis of IL-6 Citrullination. IL-6 and citrullinated IL-6 were reacted with Cayman's Citrulline-specific Probe-biotin (Item No. 17450) and detected using Streptavidin:HRP (Item No. 16747).

10	20	30	40	50
GVP PGESKD	VAAPH RQPLT	SSERIDKQIR	YLDGISALR	KETCNKSNMC
60	70	80	90	100
ESSKEALAE	NLNLPKMAEK	DGCFQSGFNE	ETCLVKIITG	LLEFEVYLEY
110	120	130	140	150
LQNRFSSEE	QARAVQMSTK	VLIQFLQKKA	KNLDAITTPD	PTTNASLLTK
160	170	180		
LQAQNQLWLD	MTTHLILRSF	KEFLQSSLR	LRQM	

Identification of modified sites in Citrullinated IL-6 (Item No. 30173). Citrullinated IL-6 was detected by LC-MS/MS and analyzed using Mascot and Scaffold PTM software. Deiminated arginines are indicated in teal.

Citrullination sites shown are representative of typical results. Batch-to-batch variations may occur.

**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, 07/12/2022

**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-6 (IL-6) is a cytokine encoded by the *IL6* gene in humans with roles in both the initiation and resolution of inflammatory responses.<sup>1,2</sup> IL-6 is produced in response to an inflammatory event, induces recruitment of neutrophils to the site of inflammation, and signals through the membrane-bound IL-6 receptor (IL-6R) to upregulate the synthesis of acute phase proteins, lipolysis in the liver, and induced differentiation of T and B cells. Following initiation of this classical signaling pathway, a soluble form of IL-6R (sIL-6R) is released from neutrophils to promote recruitment of monocytes and macrophages, induction of fever *via* IL-6 action in the hypothalamus, and production of anti-inflammatory M2 macrophages to initiate tissue repair. Production of IL-6 is dysregulated in various chronic inflammatory diseases, including rheumatoid arthritis, Castleman disease, polymyalgia rheumatica, and giant cell arteritis.<sup>3</sup> It also induces tumor growth and metastasis *via* activation of JAK/STAT signaling.<sup>4</sup> Recombinant human IL-6 can be citrullinated at arginine residues by protein arginine deiminase 4 (PAD4) *in vitro*.<sup>5</sup> Citrullinated IL-6 and anti-citrullinated peptide antibodies (ACPAs) reactive with citrullinated IL-6 have been found in synovial fluid and sera, respectively, from patients with rheumatoid arthritis. This product contains purified IL-6 (human, recombinant) (Item No. 30174) that has been modified with PAD4 enzyme, which is subsequently depleted by affinity chromatography.

## References

1. 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).
2. Rose-John, S. Interleukin-6 family cytokines. *Cold Spring Harb. Perspect. Biol.* **10**(2), (2018).
3. Schett, G. Physiological effects of modulating the interleukin-6 axis. *Rheumatology (Oxford)* **57**(suppl 2), ii43-ii50 (2018).
4. Lacina, L., Brábek, J., Král, V., *et al.* Interleukin-6: A molecule with complex biological impact in cancer. *Histol. Histopathol.* **34**(2), 125-136 (2019).
5. Lu, C., Ohara, R., Campbell, P., *et al.* Citrullination of interleukin 6 augments its pro-inflammatory capacity and signaling potency through interleukin-6 receptor in rheumatoid arthritis [abstract]. *Arthritis Rheumatol.* **71** (suppl 10) (2019).

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