

IL-8 (human), (recombinant)

Interleukin-8 (IL-8) is a chemokine, a member of the cytokine family that displays chemotactic activity for specific types of leukocytes. IL-8 regulates leukocyte recruitment and trafficking and also has mitogenic and angiogenic properties. IL-8 expression has been demonstrated in neuroendocrine tumor cells in human cancer tissue.

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Interleukin-8, CXCL8, Monocyte-derived neutrophil chemotactic factor, MDNCF, Neutrophil activating factor, NAF, NAP-1
Formulation	Lyophilized with no additives.
MW	~8.4 kDa
Purity	≥98% (SDS & HPLC)
Purity Detail	Sterile filtered through a 0.2 micron filter.
Source	Produced in <i>E. coli</i> .
UniProt ID	P10145

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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