

Human Interleukin 22 (Human Cell- Expressed)

Ordering Information

[Order Online »](#)

PBL-11701-1	10µg
PBL-11701-2	50µg

Manuals, SDS & CofA

[View Online »](#)

- Recombinant Human IL-22 protein with high purity
- Authentic human cell-expressed with no tag
- Suitable for Direct Binding ELISA, Western Blot and IL-10 Secretion Bioassay

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IL-22

Application Notes Tested Applications: Direct Binding ELISA, Western Blot, IL-10 Secretion Bioassay

Biological Activity Measured by IL-10 induction in COLO 205 human colorectal cells.

Endotoxin Content < 1 EU/μg

Formulation Phosphate buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

MW 16.8 kDa

Purity ≥ 95% (SDS-PAGE)

Purity Detail A combination of ion exchange and size exclusion chromatography.

Source Recombinant DNA expressed in Human Embryonic Kidney Cells (HEK293).



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

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