

Perforin 1 (human), (recombinant) (aa 22- 555)

Perforin 1 (PRF1) is a pore forming cytolytic protein found in the granules of cytotoxic T lymphocytes (CTLs) and NK cells. Upon degranulation, perforin binds to the target cell's plasma membrane, and oligomerizes in a Ca^{2+} dependent manner to form pores on the target cell. The pore formed allows for the passive diffusion of a family of pro-apoptotic proteases, known as the granzymes, into the target cell.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-PRT313-0010	10µg
ENZ-PRT313-0200	200µg

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE of active human recombinant PRF1

Handling & Storage

Use/Stability	Stable at 2-8°C for one month. Stable at -80°C for up to 12 months.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Lymphocyte pore forming protein, PFP, Cytolysin, PRF1
Formulation	Lyophilized from 20mM Tris, 150mM NaCl, pH 8.0, containing 0.01% skl, 5% Trehalose.
MW	62kDa
Purity	≥80%
Reconstitution	Reconstitute in ddH ₂ O to a concentration of 0.1-0.5 mg/mL. Do not vortex.
Source	Produced in <i>E. coli</i> . Recombinant human perforin 1 (PRF1) (aa 22-555).

Last modified: January 13, 2025



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-
fr@enzolifesciences.com](mailto:info-fr@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)