# 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<sup>2+</sup> 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 »

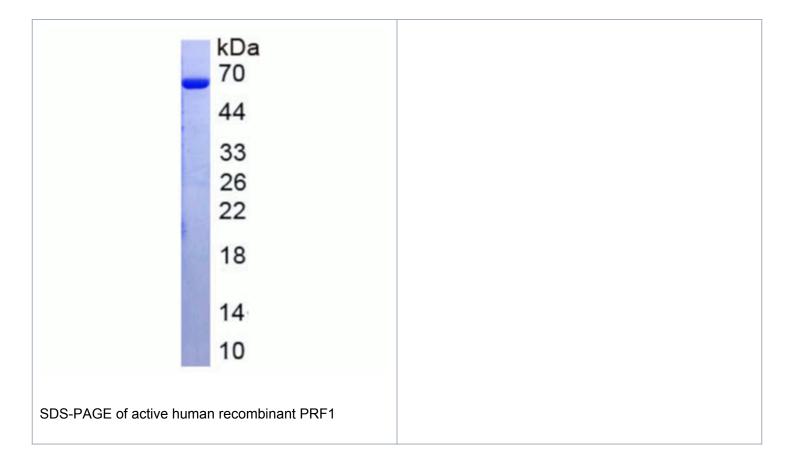
## **Ordering Information**

**Order Online** »

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

Manuals, SDS & CofA

View Online »



### **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 ddH2O to a concentration of 0.1-0.5

mg/mL. Do not vortex.

**Source** Produced in *E. coli*. Recombinant human perforin 1

(PRF1) (aa 22-555).

Last modified: January 13, 2025

