

# PRODUCT INFORMATION



## Histidine-rich Glycoprotein (human, recombinant)

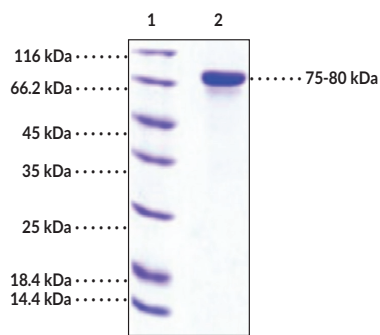
Item No. 37017

### Overview and Properties

<b>Synonyms:</b>	Histidine-proline-rich Glycoprotein, HPRG, HRG
<b>Source:</b>	Active recombinant human C-terminal His-tagged histidine-rich glycoprotein expressed in HEK293 cells
<b>Amino Acids:</b>	19-525
<b>Uniprot No.:</b>	P04196
<b>Molecular Weight:</b>	59 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥97% estimated by SDS-PAGE
<b>Supplied in:</b>	Lyophilized from sterile PBS, pH 7.4
<b>Endotoxin Testing:</b>	<1.0 EU/μg, determined by the LAL endotoxin assay
<b>Bioactivity:</b>	See figures for details

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

### Image



Lane 1: MW Markers  
Lane 2: Histidine-rich Glycoprotein

**SDS-PAGE Analysis of Histidine-rich Glycoprotein.** This protein has a calculated molecular weight of 59 kDa. It has an apparent molecular weight of approximately 75-80 kDa by SDS-PAGE under reducing conditions due to glycosylation.

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, 01/27/2023

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

---

Histidine-rich glycoprotein is an abundant plasma glycoprotein that is involved in diverse biological processes, including angiogenesis, hemostasis, cell adhesion, and immunity.<sup>1,2</sup> It is composed of a signal peptide and two cystatin-like regions in the N-terminal domain, a central domain with a histidine-rich region (HRR) and two proline-rich regions (PRRs) on either side, and a C-terminal domain. Histidine-rich glycoprotein is mainly produced by the liver but is also found in platelets and megakaryocytes.<sup>3,4</sup> It has many binding partners, including IgG, immunoglobulin  $\gamma$  receptor (Fc $\gamma$ R), and complement 1q (C1q), which bind to the N-terminal domain, zinc, heme, and microbes, which bind to the HRR, heparin and heparan sulfate, which bind to the N-terminal domain and HRR, plasminogen and thrombospondin, which bind to the N-terminal and C-terminal domains, and phospholipids, fibrinogen, and complement components, which have undefined binding regions.<sup>2</sup> Histidine-rich glycoprotein is also able to bind to various cell types such as T cells, B cells, erythrocytes, endothelial cells, and several cancer types. It modulates macrophage polarization, facilitates the clearance of apoptotic and necrotic cells, promotes antitumor immunity, and has antibacterial and antiviral activities.<sup>1,2</sup> Overexpression of *HRG* decreases tumor burden, metastasis, hypoxia, and blood vessel density and increases tumor-associated macrophage (TAM) accumulation in several mouse xenograft models.<sup>5</sup> Increased plasma levels of histidine-rich glycoprotein are associated with the development of post-thrombotic syndrome following a venous thromboembolic event (VTE).<sup>6</sup> Cayman's Histidine-rich Glycoprotein (human, recombinant) protein can be used for cell-based adhesion assays. This protein consists of 518 amino acids, has a calculated molecular weight of 59 kDa, and a predicted N-terminus of Val19 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 75-80 kDa due to glycosylation.

## References

---

1. Pan, Y., Deng, L., Wang, H., *et al.* Histidine-rich glycoprotein (HRGP): Pleiotropic and paradoxical effects on macrophage, tumor microenvironment, angiogenesis, and other physiological and pathological processes. *Genes Dis.* **9(2)**, 381-392 (2022).
2. Poon, I.K.H., Patel, K.K., Davis, D.S., *et al.* Histidine-rich glycoprotein: the Swiss Army knife of mammalian plasma. *Blood* **117(7)**, 2093-2101 (2011).
3. Koide, T., Foster, D., Yoshitake, S., *et al.* Amino acid sequence of human histidine-rich glycoprotein derived from the nucleotide sequence of its cDNA. *Biochemistry* **25(8)**, 2220-2225 (1986).
4. Leung, L.L.K., Harpel, P.C., Nachman, R.L., *et al.* Histidine-rich glycoprotein is present in human platelets and is released following thrombin stimulation. *Blood* **62(5)**, 1016-1021 (1983).
5. Rolny, C., Mazzone, M., Tugues, S., *et al.* HRG inhibits tumor growth and metastasis by inducing macrophage polarization and vessel normalization through downregulation of PlGF. *Cancer Cell* **19(1)**, 31-44 (2011).
6. Siudut, J., Natorska, J., Son, M., *et al.* Increased levels of histidine-rich glycoprotein are associated with the development of post-thrombotic syndrome. *Sci. Rep.* **10(1)**, 14419 (2020).

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