

Data Sheet

GSTP1, His-FLAG-Tags

Human, Recombinant, N-terminal His-FLAG-tags

Catalog #: 79415

Lot #: 190805-1

Conc.: 0.55 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 10 mM Imidazole and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

Description: Human Glutathione S-Transferase Pi 1 (GSTP1), also known as GST Class-Pi, GST3, Fatty Acid Ethyl Ester Synthase III and Deafness X-Linked 7 (DNF7), Genbank Accession No. NM_000852, a.a. 2-211(end) with N-terminal His-FLAG-tags, expressed in an *E. coli* expression system. MW=26 kDa.

References:

1. Cairns, P., *et al.*, *Clinical Cancer Research.*, 2001 Sep; **7(9)**:2727-2730.
2. Lin, X., *et al.*, *Am. J. of Pathology.* 2001 Nov; **159(5)**:1815-1826.

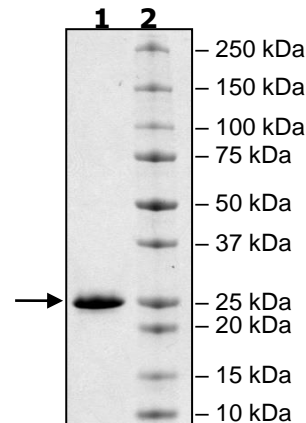
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg GSTP1
Lane 2:
Protein Marker

MW: 26 kDa

Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com