

GAS6 (human), (recombinant) (His- tag)

The growth arrest-specific 6 gene (GAS6) is a vitamin K-dependent protein that plays a role in the survival, proliferation, migration, differentiation, adhesion, and apoptosis of cells. The growth arrest-specific 6 (GAS6) has been implicated in systemic inflammation, coagulation, and plays a role in tumor progression by regulating growth in many cancers. GAS6, expressed by osteoblasts in the bone marrow, plays a significant role in the regulation of PCa cell survival during chemotherapy, which will have important implications for targeting metastatic disease. The GAS6/TYRO3-AXL-MERTK (TAM) signaling pathway is essential for full and sustained platelet activation, as well as thrombus stabilization. Inhibition of this pathway decreases platelet aggregation, shape change, clot retraction, aggregate formation under flow conditions, and surface expression of activation markers. GAS6 signaling regulates invasion, proliferation, chemotherapy-induced apoptosis of prostate cancer (PCa) cells, and GAS6 secreted from osteoblasts in the bone marrow environment plays a critical role in establishing prostate tumor cell dormancy.

Ordering Information

[Order Online »](#)

ENZ-PRT326-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

kDa

M

116

66.2

45.0

35.0

25.0

18.4

14.4



It migrates as approximately 43.9 kDa band in SDS-PAGE under reducing conditions.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Growth arrest-specific protein 6, AXL receptor tyrosine kinase ligand

Endotoxin Content <1 EU/μg (LAL method)

Formulation Lyophilized from sterile 20 mM Tris, 150 mM NaCl, 10% Glycerol, 10% Trehalose, 0.1 M Arg, pH 8.0.

MW ~43.9 kDa (SDS-PAGE)

Purity ≥90% (SDS-PAGE)

Reconstitution Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

Source Produced in HEK293 cells. A DNA sequence encoding the human GAS6 (Asp279-Ala678) was expressed with a polyhistidine tag at the N-terminus.

UniProt ID Q14393-2



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)