
STAT3 Protein

(Catalog No. E62006)

Description

STAT3 Human Recombinant encoding aa 50-240 with GST tag at N-terminal.

Molecular Weight (kDa)

49 kDa

Formulation

50mM Tris-Acetate, pH7.5, 1mM EDTA, 20% Glycerol

Source

Expressed in E.coli expression system

Size

10 µg

Storage Temp & Conditions

1 week at 14°C. For longer periods of time store vial at -20°C to -80°C. When stored at the recommended temperature, this protein is stable for 12 months. Avoid multiple freeze-thaw cycles.

Purification Notes

SDS-PAGE Purified

Application Notes

ELISA, Inhibition Assay, WB

Ordering Information:

Products	Size	Cat. No.
STAT3 Protein	10 µg	E62006 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.