

WASP-121^{GST}

(Wiskott-Aldrich Syndrome Protein)

Human, Recombinant, *E. coli*

Cat. No.	Amount
PR-369	0.25 mg

Liquid. Supplied as a 1 mg/ml solution in 50 mM Tris-HCl, pH 7.5, 100 mM NaCl, and 2 mM DTE.

The Wiskott-Aldrich syndrome protein (WASP) is the founding member of a family of proteins that have emerged as crucial effectors of Rho GTPases and activators of the cytoskeletal organizing complex Arp2/3. WASP has been shown to be intimately involved in many pathways that influence the function of the immune system. WASP-121 comprises the amino acids 201-321 of the native WASP sequence. This domain is essential for the binding of Cdc42 to WASP.

The GST-Tag facilitates the protein's application in typical GST pull down assays.

AVOID FREEZE/THAW CYCLES!

For in vitro use only!

Purity: > 95% by SDS-PAGE

Store: -80 °C

Selected References:

Thrasher A.J. (2002) WASp in immun-system organization and function. *Nature Reviews Immunology* **2**:635.

Rudolph *et al.* (1998) The Cdc42/Rac interactive binding region of the Wiskott-Aldrich-Syndrome Protein (WASP) is necessary but not sufficient for tight binding to Cdc42 and structure formation. *J. Biol. Chem.* **273**:18067.