

## BUB1 Protein

Recombinant protein expressed in Sf9 cells

**Catalog # B11-35G**

Lot # S179-4

### Product Description

Recombinant human BUB1(756-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_004336](#).

### Gene Aliases

BUB1A, BUB1L, hBUB1

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

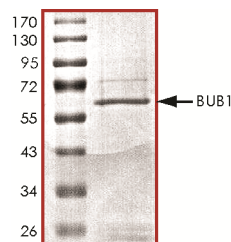
### Scientific Background

BUB1 is a kinase which involved in spindle checkpoint function that in part by phosphorylating a member of the mitotic checkpoint complex and activating the spindle checkpoint. BUB1 is highly expressed in thymus, ovary, and testis (1). BUB1 maintains the steady-state level and centromeric localization of SGO1, and they concluded that BUB1 protects centromeric cohesion in mitosis through SGO1 (2). BUB1-mediated phosphorylation of CDC20 is required for proper checkpoint signaling and that inhibition of APC by BUB1 may partly account for the sensitivity of the spindle checkpoint. BUB1 kinase creates a mark for shugoshin localization and the correct partitioning of chromosomes.

### References

1. Leland, S.et.al: Heterozygosity for a Bub1 mutation causes female-specific germ cell aneuploidy in mice. Proc. Nat. Acad. Sci. 106: 12776-12781, 2009.
2. Tang, Z.et.al: Human Bub1 protects centromeric sister-chromatid cohesion through Shugoshin during mitosis. Proc. Nat. Acad. Sci. 101: 18012-18017, 2004.

### Purity



The purity of BUB1 was determined to be **>80%** by densitometry. Approx. MW **63 kDa**.

## BUB1 Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	B11-35G
Specific Lot Number	S179-4
Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**