

## FOXL2 Protein

Full-length human recombinant protein expressed in Sf9 cells

**Catalog # F71-30BG**

Lot # B2024-8

### Product Description

Full-length recombinant human FOXL2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_023067](#).

### Gene Aliases

BPES; BPES1; PFRK; PINTO; POF3

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . 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

Fork-head transcriptional factor forkhead box L2 (FOXL2) is a member of the large family of Forkhead Box (Fox) domain transcription factors. Plays a key role in sex differentiation and oogenesis. FOXL2 is involved in virtually all stages of ovarian development and function, as well as in granulosa cell (GC)-related pathologies. It maintains GC identity through the repression of testis-specific genes.

### References

1. Baron D, et al: Foxl2 gene and the development of the ovary: a story about goat, mouse, fish and woman. *Reprod Nutr Dev* 45:377-382 (2005).
2. Georges A, et al: FOXL2: a central transcription factor of the ovary. *J Mol Endocrinol*. 2013 52(1):R17-33.

### Purity



The purity of FOXL2 was determined to be **>70%** by densitometry. Approx. MW **~73kDa**.

## FOXL2 Protein

Full-length human recombinant protein expressed in Sf9 cells

Catalog #	F71-30BG
Lot #	B2024-8
Purity	>70%
Concentration	0.05 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . 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.**