

## DPY30 Protein

Full-length recombinant protein expressed in E. coli cells

**Catalog # D369-30H**

Lot # B2178-6

### Product Description

Recombinant full-length human DPY30 was expressed in E.coli cells using an N-terminal His tag. The gene accession number is [NM\\_032574](#).

### Gene Aliases

Cps25; HDPY-30; Saf19

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

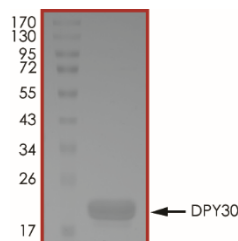
### Scientific Background

As a core subunit of the SET1/MLL histone methyltransferase complexes, DPY30 modulates H3K4 methylation in vitro, and directly regulates chromosomal H3K4 trimethylation (H3K4me3) throughout the mammalian genome. The methylation event has been shown to affect embryonic stem cells (ESCs) maintenance and differentiation. Depletion of Dpy-30 does not affect ESC self-renewal, but significantly alters the differentiation potential of ESCs, particularly along the neural lineage.

### References

1. Jiang H, et al: Role for Dpy-30 in ES cell-fate specification by regulation of H3K4 methylation within bivalent domains. Cell. 2011 Feb 18;144(4):513-25.
2. Patel A, et al: On the mechanism of multiple lysine methylation by the human mixed lineage leukemia protein-1 (MLL1) core complex. J Biol Chem. 2009 Sep 4;284(36):24242-56.

### Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **20kDa**.

## DPY30 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog #	D369-30H
Lot #	B2178-6
Purity	>95%
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.**