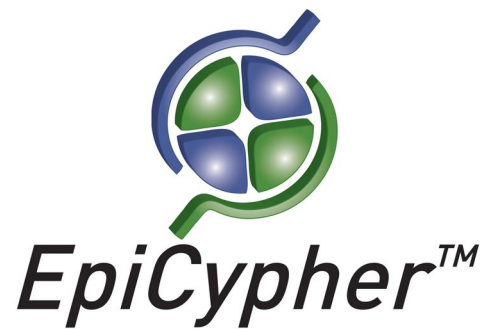


# RAG2 PHD Domain, Recombinant

**Catalog No.** 15-0066  
**Lot No.** 17025001  
**Pack Size** 100 µg



**Type** PHD                      **Expressed In** *E. coli*  
**Mol. Wgt.** 43 kDa              **Epitope Tag** GST

## Product Description:

RAG2 PHD Domain, Recombinant (V(D)J Recombination Activating Protein 2, accession number Q12830, amino acids 387 - 493) containing an N-terminal GST tag, expressed in *E. coli*. RAG2 is part of a protein complex that mediates DNA cleavage during immunoglobulin gene V(D)J recombination. RAG2 contains a PHD domain that interacts with H3K4me3, which recruits the RAG complex to sites of active chromatin.

## Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 50 mM HEPES pH 7.3, 200 mM NaCl, 1 mM DTT, 20% glycerol.

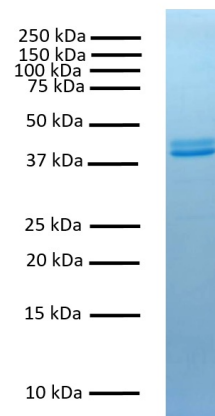
## Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

## Application Notes:

RAG2 PHD Domain, Recombinant is useful for protein binding and screening experiments examining methylated protein and peptide substrates.

## References:



**Protein Gel Data:** RAG2 PHD Domain, Recombinant Human (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

This product is for *in vitro* research use only and is not intended for use in humans or animals.