

SAP145, GST-Tag

Human, recombinant, N-terminal GST-tag

Catalog #: 79111

Lot #: 160817-1

Conc.: 0.71 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04%
Tween20, 16mM glutathione, 20% glycerol

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles.*

References:

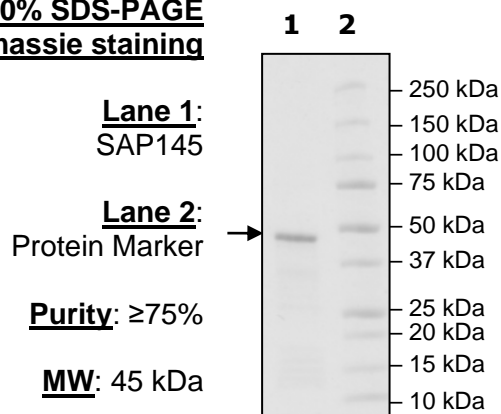
1. Yang, Y., *et al.* (2015) *Nat Commun* **6**, 6428
2. Falk, S., *et al.* (2016) *Nat Commun* **7**, 13573

Description: Human SAP145, a substrate for PRMT9, also known as Splicing Factor 3b Subunit 2 (SF3B2), GenBank accession No. NM_006842, a.a. 401-550 with an N-terminal GST-tag, expressed in an *E. coli* expression system. MW = 45 kDa

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com