

Data Sheet

Histone H3, His-Tag

Human, recombinant, N-terminal His-tag

Catalog #: 52023

Lot#: 131122 **Conc.:** 1.19 mg/ml

Formulated in: 8 mM PBS pH 7.4, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 20% glycerol.

Stability: >6 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Loyola, A. *et al.*, *Mol. Cell* **24** (2), 309-316 (2006).
2. Hans, F. and Dimitrov, S. *Oncogene* **20**(24), 3021-3027 (2001).

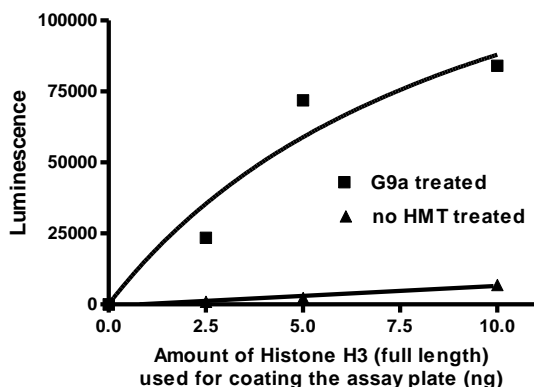
Description: Human Histone 3 (HIST1H3E), GenBank Accession No. NM_003532, a.a. 2-136(end) with N-terminal His-tag MW = 15.4 kDa, expressed in an *E. coli* expression system.

Assay Conditions: Histone H3 was coated onto a 96-well plate, and methylation by G9a was carried out at room temperature for 60 min. Methylated Histone H3 was detected using a chemiluminescence assay.

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

Quality Assurance

Histone H3, Methylation Assay

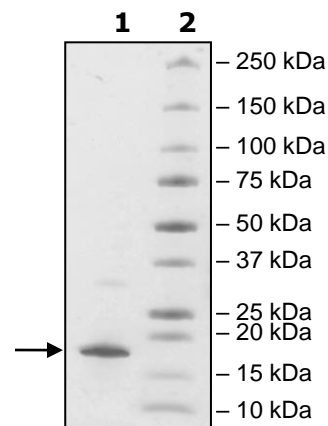


4-20% SDS-PAGE Coomassie staining

Lane 1:
2 µg Histone H3

Lane 2:
Protein Marker

MW: 15 kDa
Purity: ≥90%



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