

## Data Sheet

### $\beta$ -glucosyltransferase

T4 phage, recombinant, C-terminal His-tag

**Catalog #:** 72000

**Lot #:** 110111      **Conc:** 1.66mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, 20% glycerol.

**Stability:** >6 months at  $-80^{\circ}\text{C}$

#### References:

1. Song, C.X., *et al.* (2011). *Nat Biotechnol.* **29(1)**: 68-72.
2. Kornberg, S.R., *et al.* (1961). *J. Biol. Chem.* **236**: 1487-1493.

#### Description:

T4-phage beta-glucosyltransferase, also known as UDP glucose-DNA  $\beta$ -glucosyltransferase, Genbank Accession No. NP\_049658, a.a. 1-351(end) with C-terminal His-tag, MW= 41.6 kDa, expressed in *E. coli*.

#### Application:

Useful for the differentiation of hydroxymethylcytosine (HMC) from methylcytosine in DNA, via glucosylating HMC and protecting HMC from endonuclease cleavage.

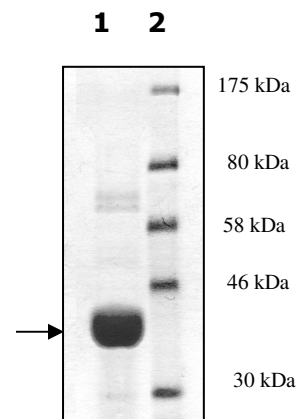
### Quality Assurance

#### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
8  $\mu\text{g}$   $\beta$ -glucosyltransferase

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 42 kDa  
**Purity:**  $\geq 83\%$



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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)