

# **EP300** Protein

(Catalog No. E24000)

### Description

Human E1A binding protein p300, also known as EP300, KAT3B, or RSTS2, GenBank Accession No. NM\_001429, a.a. 965-1810, with N-terminal His- GST-tag.

#### Molecular Weight (kDa) 125 kDa

125 KDa

#### Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, and 20% glycerol.

#### Source

Expressed in Sf9 cells via a baculovirus expression system.

## Size

50 µg

Storage Temp & Conditions >6 months at -80°C

>6 months at  $-80^{\circ}$ C

**Purification Notes** Glutathione Sepharose

#### **Application Notes**

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

#### Ordering Information:

Products	Size	Cat. No.
EP300 Protein	50 µg	E24000 -1