

Product Information

Description:	Recombinant human BRDT (bromodomain testis-specific) protein, encompassing amino acids 22-382, corresponding to the bromodomain (BD) 1 and 2. This construct contains an N-terminal GST-Tag followed by a PreScission protease cleavage site. The recombinant protein was affinity purified.
Background:	BRDT (bromodomain testis-specific) protein, also known as CT9 and BRD6 (bromodomain-containing protein 6), is a key regulator of gene expression and chromatin organization. This protein is a testis-specific member of the bromodomain and extra-terminal motif (BET) family and is important for male meiosis and spermatogenesis. BRDT possesses two bromodomain motifs (BD1 and BD2) at its N-terminus and a C-terminal extra-terminal (ET) domain. Dysregulation of this protein is associated with male infertility and several types of cancer including lung, ovarian, and esophageal squamous cell carcinoma. BET bromodomain inhibitors, such as JQ1, are under development as potential cancer therapeutics.
Species:	Human
Construct:	BRDT (BD1+BD2) (GST-Pre-22-382)
Concentration:	0.57 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.
MW:	68.1 kDa
Genbank Accession:	NM_207189
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful for SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

