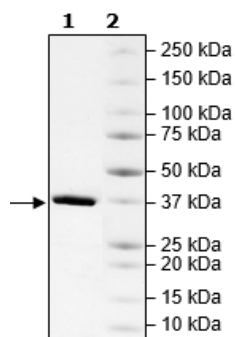


Product Information

Description:	Recombinant human bromodomain CECR2 (cat eye syndrome chromosome region, candidate 2), encompassing amino acids 430-543, which correspond to the single bromodomain. This construct contains an N-terminal GST-tag. This protein was affinity purified.
Background:	CECR2 (cat eye syndrome chromosome region candidate 2) is an epigenetic factor containing a bromodomain able to recognize acetylated lysines. It is involved in DNA Damage Responses (DDR), neurulation, spermatogenesis, and reprogramming. It has been linked to cancer and was identified as an epigenetic driver of metastatic breast cancer. It can form a complex with RELA (v-rel avian reticuloendotheliosis viral oncogene homolog A), increasing the access to chromatin of EMT (epithelial-mesenchymal transition) related genes and NF- κ B (nuclear factor kappa B) target genes. The use of CECR2 inhibitors, such as GNE-886 and NVS-CECR2-1, is a promising area in cancer therapy.
Species:	Human
Construct:	CECR2 (GST-430-543)
Concentration:	3.70 mg/ml
Expression System:	<i>E. coli</i>
Purity:	$\geq 90\%$
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris, pH 8, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween, and 20% glycerol
MW:	40 kDa
Genbank Accession:	NM_031413
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.
Assay Conditions:	Assay was performed according to CECR2 TR-FRET Assay Kit (BPS Bioscience #32622) with various amounts of CECR2.
Applications:	Useful for binding studies.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CECR2 Binding

