

Data Sheet

CECR2

Human, recombinant, N-terminal His tag

Catalog #: 31046

Lot #: 141009-G **Conc.:** 0.44 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl. 0.04% Tween-20 and, 20% glycerol

Stability: >6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References

1. Chen J, et al., *Int J Dev Biol.* 2010; **54(5)**: 925-9.
2. Lee SK, et al., *Mol Cells.* 2012; **34(1)**: 85-91.

Description: Human Bromodomain CECR2 (cat eye syndrome chromosome region, candidate 2), amino-acids 430-543 (GenBank Accession No. NM_031413) with N-terminal His-tag, MW = 14.4 kDa, expressed in an *E. coli* expression system.

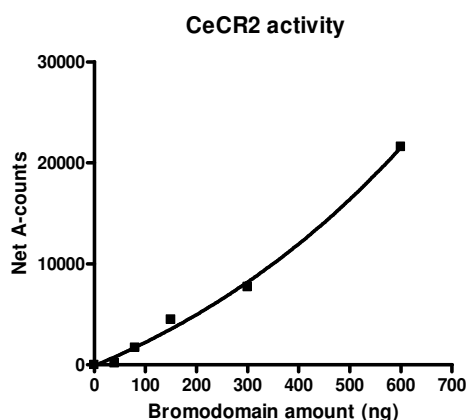
Assay Conditions: CECR2 was incubated with various amounts of BET ligand in assay buffer in a 10- μl reaction for 30 min at RT. Reactions with H4 1-21[B] were prepared in parallel as control. AlphaLISA® Ni acceptor beads (1:250) were added and incubation continued for 30 min more. Finally, Streptavidin Donor beads (1:125) were added and mixture was incubated for 10 min at RT before A-counts were measured.

Application:

Useful for the study of bromodomain binding assays, screening inhibitors and selectivity profiling.

Quality Assurance

Specific Activity

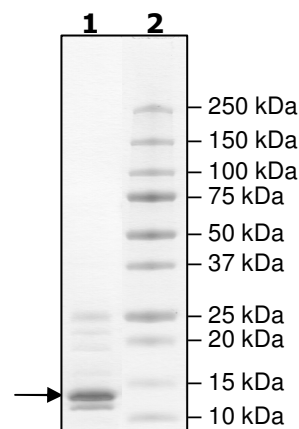


4-20% SDS-PAGE Coomassie staining

Lane 1:
1.9 μg CECR2

Lane 2:
Protein Marker

MW: 14.4 kDa
Purity: $\geq 70\%$



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

**MHHHHHHTKDLFELDDFTAMYKVL DVVKAHKDSWPFLEPVDESYAPNYYQIIKAPM
DISSMEKKLNGGLYCTKEEFVNDMKT MFRNCRKYNGESSEYTKMSDNLERC FHRAM
MKHFPGED**

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