

CRE/CREB – GLuc-HRP - cAMP Reporter

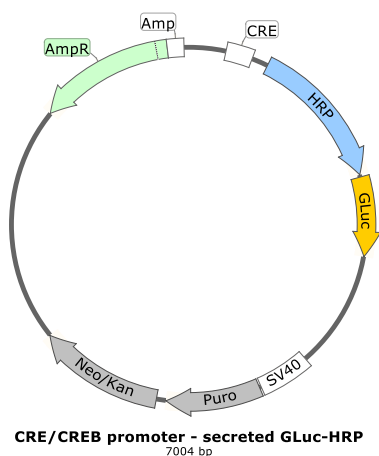
by InTouch BioSolutions



This cAMP reporter contains the control of multimerized cAMP response element (CRE) located upstream of the secreted Gaussia Luciferase (wt) - HRP (horseradish peroxidase) gene fusion. Expression of the GLuc-HRP fusion is induced by intracellular cAMP as a result of binding of the cAMP response element binding protein (CREB) to CRE. Secreted GLuc-HRP can be measured in the supernatant (see data below).

Plasmid details

- Response element: cAMP (cyclic adenosine monophosphate)
- Promoter containing CRE sites for CREB binding
- Reporter: Secreted Gaussia Luciferase (wt) - HRP (horseradish peroxidase) fusion protein
- Selection markers: SV40 promoter – Puromycin, Neomycin and Kanamycin resistance genes
- Plasmid size: ~7 kb (contains bacterial-derived sequences)
- Storage: -20°C. For long-term storage, the recommended storage is -80°C.



Application

- Expressed Gaussia Luciferase – HRP is secreted into the supernatant, thus cell lysis is not necessary to measure activity. Ideal for multiple sampling in time-dependent assays.
- Transcription regulation, virus-cell interactions, compound screening, post-translational modifications, GPCR signaling, cell signaling, promoter analysis.
- End-user preference on bioluminescence or colorimetric detection

Sample protocol

Approximately 1 million human melanoma cells were transfected with 10ug of CRE/CREB – GLuc-HRP plasmid by calcium phosphate method and seeded in a 96-well (5000 cells/well) flat bottom plate. The next day, cells were incubated with a GPCR agonist (alpha-MSH; 5nM) mixed with increasing amounts of inhibitor.

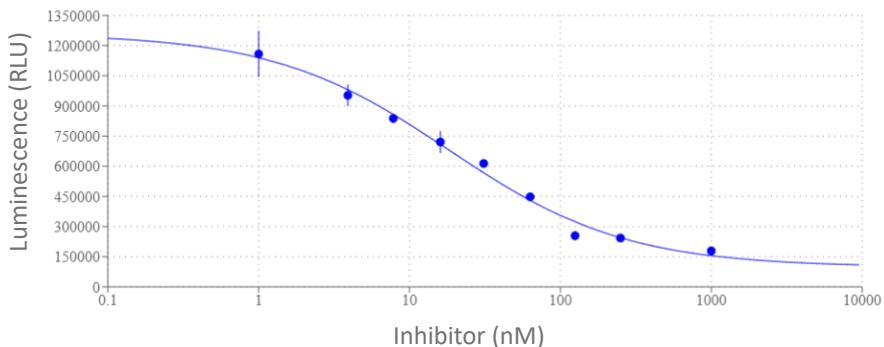
CRE/CREB – GLuc-HRP - cAMP Reporter

by InTouch BioSolutions



Data

A. Gaussia Luciferase activity



IC_{50} (nM) = 18.4

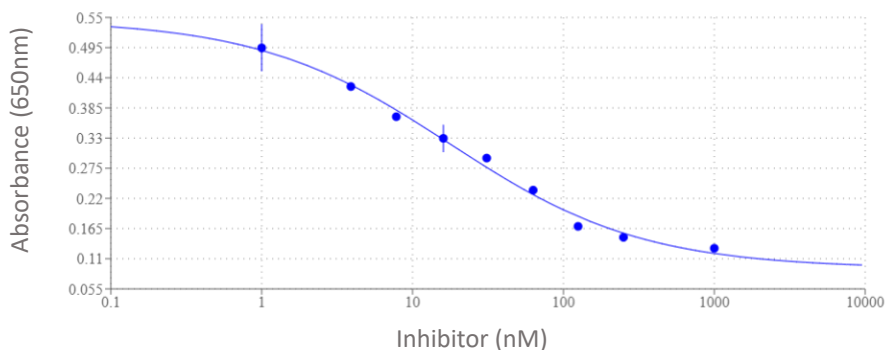
Max: 1,272,648

Min: 177,490

Error bar: SEM from duplicate wells

Figure A) After 24 hours, 10 μ l from each well was added to wells containing native CTZ (Cat. #303, Nanolight Technology) and luminescence (RLU) measured using Molecular Devices Spectramax L plate reader.

B. HRP activity



IC_{50} (nM) = 17.7

Max: 0.537

Min: 0.127

Error bar: SEM from duplicate wells

Figure B) After 48 hours, 5 μ l from each well was added to wells containing TMB substrate and the absorbance measured at 650 nm using Molecular Devices Thermomax plate reader. Inset is a representative image from the duplicate wells showing

Terms

This product is intended to be used only for research use and is not intended for diagnostic or therapeutic use or administration to animals or humans. Please send an email to info@intouchbio.com to obtain the material transfer agreement (MTA) for further details regarding the use of this product.