

CRHR1 Polyclonal Antibody

(Catalog # A53655)

Background

G-protein coupled receptor for CRH (corticotropin-releasing factor) and UCN (urocortin). Has high affinity for CRH and UCN. Ligand binding causes a conformation change that triggers signaling via guanine nucleotide-binding proteins (G proteins) and down-stream effectors, such as adenylate cyclase. Promotes the activation of adenylate cyclase, leading to increased intracellular cAMP levels. Inhibits the activity of the calcium channel CACNA1H. Required for normal embryonic development of the adrenal gland and for normal hormonal responses to stress. Plays a role in the response to anxiogenic stimuli.

Description

CRHR1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P34998

Purification

Protein G purified

Immunogen

Recombinant Human Corticotropin-releasing factor receptor 1 protein (24-121AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Corticotropin-releasing hormone receptor 1 CRHR1 CRFR, CRFR1, CRHR

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western blot

All lanes: CRHR1 Polyclonal Antibody at 16µg/ml

Lane 1: Hela whole cell lysate

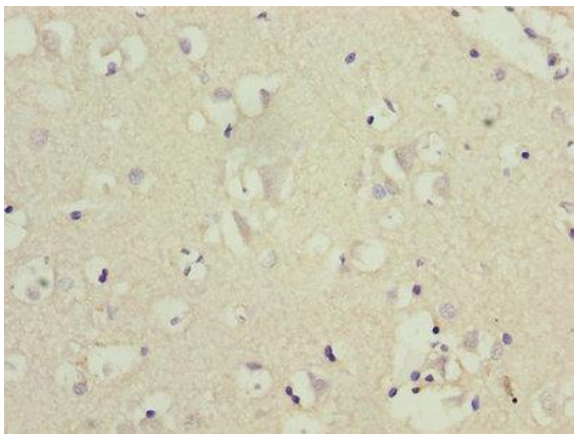
Lane 2: HL60 whole cell lysate

Secondary

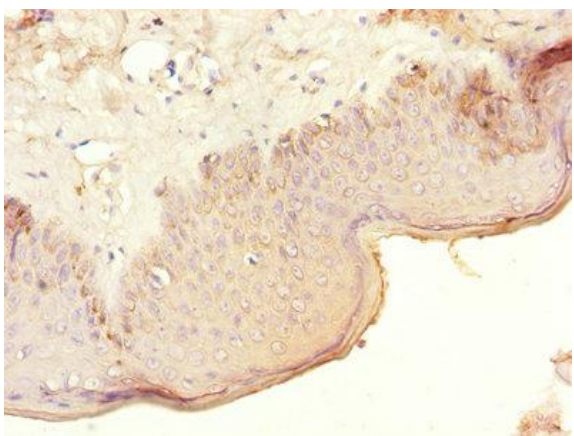
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 51, 48, 44, 47, 29 kDa

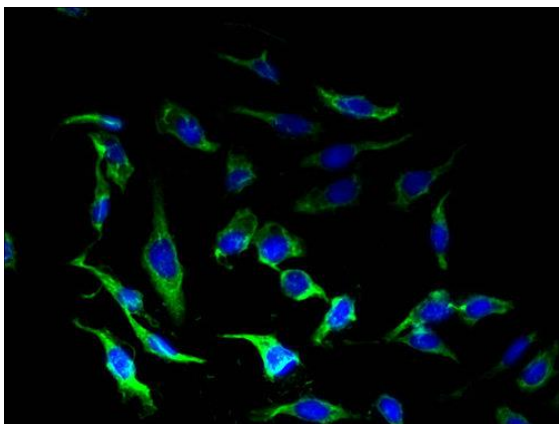
Observed band size: 51 kDa



Immunohistochemistry of paraffin-embedded human brain tissue at dilution of 1:100



Immunohistochemistry of paraffin-embedded human skin tissue at dilution of 1:100



Immunofluorescent analysis of Hela cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)