Prolactin (human), (recombinant)

Prolactin is a hormone, produced by the anterior pituitary gland, that has been shown to induce lactation in the mammary glands after birth. It has since been shown to have many diverse regulatory functions in the body including reproductive, metabolic, osmotic, immunological, and behavioral roles. Elevated levels of prolactin can be associated with hypothalamic disease, cancer, hypothyroidism, and kidney disease. Prolactin levels can be controlled inhibitory factors, such as dopamine. Prolactin has also been shown to stimulate proliferation of Nb2-11cells in a dose dependent manner.

Ordering Information

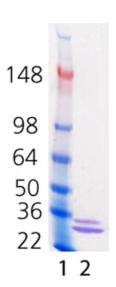
Order Online »

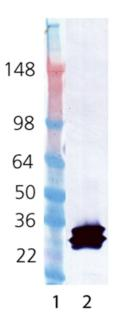
ENZ-PRT194-0050

50µg

Manuals, SDS & CofA

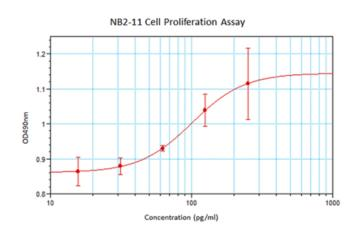
View Online »





Coomassie stained SDS-PAGE. Lane 1, Molecular weight marker. Lane 2, $1.0\mu g$ prolactin (ENZ-PRT194-0050).

Western blot analysis of prolactin probed with rabbit anti FLAG antibody: Lane 1: MWM, Lane 2: 100ng prolactin (ENZ-PRT194-0010) probed FLAG antibody



Nb2-11 cell proliferation assay after treatment with prolactin (ENZ-PRT194-0050): Nb2-11 cells were placed in stationary media at 50,000cells/well in 96 well plates. They were treated with a titered concentration of prolactin. The cells were grown at 37OC / 5% CO2 for 48 hours. Proliferation was measured using an MTT cell viability assay. The ED50 range is 30 – 200pg/ml.

Handling & Storage

Use/Stability Stable for at least 6 months after receipt when stored at -20°C.

Handling After reconstitution, prepare aliquots and store at -80°C. Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Luteotropic hormone, Luteotropin

Biological Activity Positive for NB2-11 cell proliferation.

Endotoxin Content ≤100EU/mg

Formulation Lyophilized from 1X PBS, pH 7.5.

MW Expected 24.6kDa, actual is doublet at ~35kDa and

~30kDa.

Purity ≥90% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Source Produced in HEK 293 cells. Prolactin (aa 29-227) is fused

at the N-terminus to a FLAG-tag, 13 amino acid spacer,

and a His-tag.

UniProt ID P01236

Last modified: May 29, 2024

