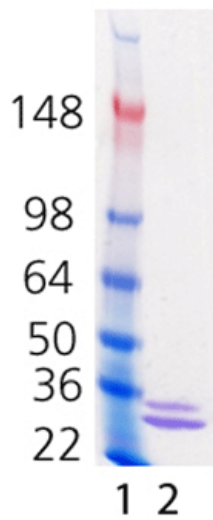


Prolactin (human), (recombinant)

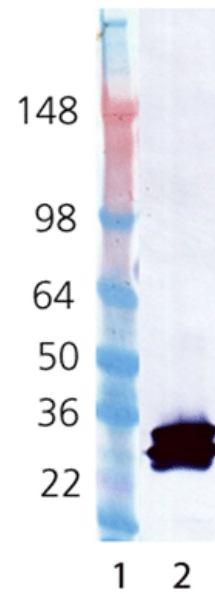
Prolactin is a hormone, produced by the anterior pituitary gland, that has been shown to induce lactation in the mammary glands after birth. It has since been shown to have many diverse regulatory functions in the body including reproductive, metabolic, osmotic, immunological, and behavioral roles. Elevated 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 Nb2-11 cells in a dose dependent manner.

Manuals, SDS & CofA

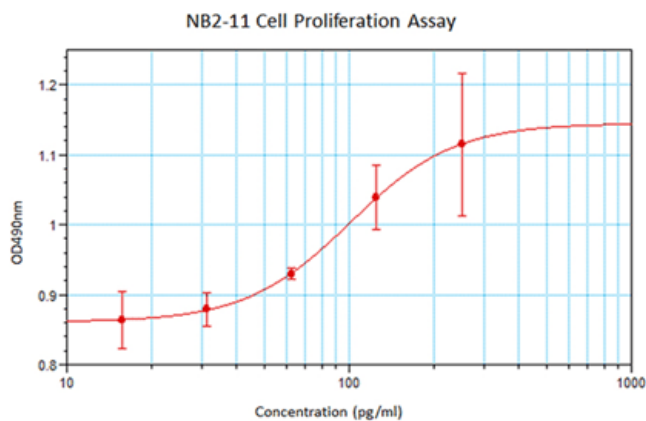
[View Online »](#)



Coomassie stained SDS-PAGE. Lane 1, Molecular weight marker. Lane 2, 1.0µg 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 titrated 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com