



6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Human Follicle Stimulating Hormone Recombinant

Catalog #: 90138
Lot #: 240207

Description: Recombinant human Follicle Stimulating Hormone, or FSH, is a heterodimeric protein consisting of 92 amino acids alpha chain and 111 amino acids beta chain. FSH migrates as an approximately 27 kDa glycoprotein.

Source: Optimized DNA sequence encoding human FSH mature chain was expressed in *CHO*.

Formulation: Lyophilized from a 0.2 µm filtered solution.

Reconstitution: A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/ml. This solution can then be diluted into other buffers.

Storage: The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots

at 2 - 8° C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

Purity: >95%, as determined by SDS-PAGE and HPLC.

Endotoxin Level: Endotoxin level was found to be < 0.1 ng/µg (1EU/µg), using the LAL gel clot method.

Biological Activity: Activity was measured to be 1.36 x 10⁴ IU/mg protein.

References:

1. *J. Endocrinol.*, Oct 2009; **203**: 167 - 174.
2. *Endocr. Relat. Cancer*, Sep 2009; **16**: 1005 – 1016.