

Certificate of Analysis

Human FGF-4, human cell-expressed (HCE)

Catalog No: 11632-2

Lot No: H00006

Expiration: One year from date of receipt

Size: 10 g

Description: HumanKine® Recombinant Human FGF-4

EC50: Typically 0.25 to 1.25 ng/ml

Formulation: Lyophilized in PBS

Endotoxin: < 1 EU/μg

Molecular Weight: 17 and 27 kDa, monomer, glycosylated

Purity: > 95%

Purity Confirmation: SDS-PAGE stained with Coomassie blue

Source: Human Cell Expressed (HCE) in HEK-293 cells in the absence of animal-derived products

Reconstitution: Briefly centrifuge the vial before opening. It is recommended to reconstitute the protein in sterile PBS containing 0.1% recombinant human serum albumin, low endotoxin (please see product catalog #11600 HSA).

Assay Used to Measure Bioactivity: FGF-4 specific activity was determined by the dose-dependent stimulation of the proliferation of the BALB/c 3T3 cell line.

References:

1. Kosaka, N., *et al* (2009) *Dev Dyn* 238, 265-276.
2. Lee, S. Y., *et al* (2000) *Molecules and Cells* 10, 684-691.

Product Information:

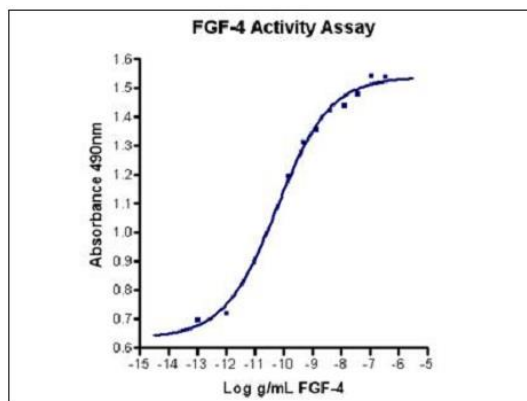


Figure 1. Human FGF-4 Activity Assay

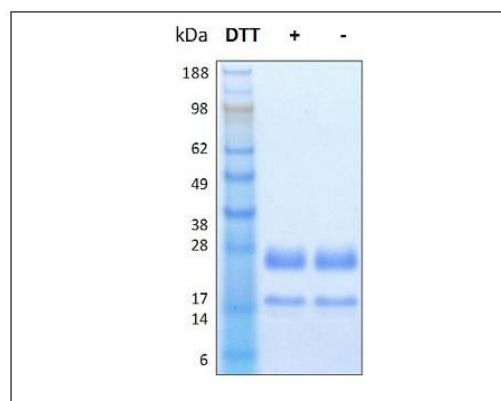


Figure 2. Purity Confirmation on SDS-PAGE



pbl assay science

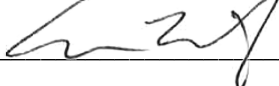
131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: info@pbl assaysci.com | W: pbl assaysci.com

Shipping Conditions: Room Temperature, 22-25°C

Physical State of Product During Shipping: Lyophilized

Storage Conditions/Comments: Lyophilized products are stable for one year at -20°C to -80°C from date of receipt. Upon reconstitution, the cytokine can be stored at 4°C for short term only (< 1 week), or at -20°C to -80°C in aliquots for long term. Avoid repeated freeze-thaw cycles. Refreezing should be performed on dry ice or in a dry ice/alcohol bath. For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____ 

Date: June 19, 2014

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.

Rev. 00