



Certificate of Analysis

Human Interleukin-22, human cell-expressed (HCE)

Catalog No: 11701-2

Lot No: 7223

Size: 50 µg/vial

Description: Recombinant Human Interleukin-22, human cell-expressed

Volume: 0.1 ml

Concentration: 0.500 mg/ml

Activity (ED₅₀): 0.45 ng/ml, measured by its ability to induce IL-10 secretion in COLO 205 human colorectal adenocarcinoma cells (Nagalakshmi, M.L. *et al.* (2004) International Immunopharmacology 4:679)

Buffer: Phosphate buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

Endotoxin: < 1 EU/µg

Molecular Weight: 16.8 kDa; due to varying degrees of glycosylation, migrates at ~25-30 kDa, as observed by reducing SDS-PAGE

Purity: > 95% as determined by SDS-PAGE

Purification Method: A combination of ion exchange, hydrophobic interaction and size exclusion chromatography

Source: Recombinant DNA expressed in Human Embryonic Kidney Cells (HEK293)

Accession #: NP_065386.1

Synonyms: TIFa; IL-22; ILTIF; IL-TIF; IL-D110; zcyto18; TIFIL-23

Tested Applications: Direct Binding ELISA, Western Blot, IL-10 Secretion Bioassay

Optimal dilutions should be determined by each laboratory for each application.

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, the product may be stored at -20°C for short-term use (≤ 6 months). For long-term storage, we recommend storing the product at -70°C or below for retention of full activity. Thaw product vial by incubation in cold tap water until just thawed – the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Refreezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA). For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____

Date: June 25, 2020

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.