



## Certificate of Analysis

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

**Catalog No:** 11701-1

**Lot No:** 7220

**Size:** 10 µg/vial

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

**Volume:** 0.1 ml

**Concentration:** 0.100 mg/ml

**Activity (ED<sub>50</sub>):** 0.80 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 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](http://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.