



## Certificate of Analysis

### Human IL-15 / sIL-15R $\alpha$ heterodimer, human cell-expressed (HCE)

**Catalog No:** 11702-2

**Lot No:** 7262

**Size:** 50  $\mu$ g/vial

**Description:** Recombinant Human IL-15 / soluble IL-15R $\alpha$  heterodimer, human cell-expressed

**Volume:** 0.1 ml

**Concentration:** 0.5 mg/ml

**Biological Activity (ED<sub>50</sub>):** <1 ng/ml (32.1 pM). Typically 0.2-1 ng/ml, measured by its dose dependent stimulation of the proliferation of CTLL-2 cells

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

**Endotoxin:** < 1 EU/ $\mu$ g

**Molecular Weight:** 31 kDa predicted (12.8 kDa for IL-15 + 18.4 kDa for sIL-15R $\alpha$ ). Heterodimer exhibits multiple bands on reducing SDS-PAGE due to varying degrees of glycosylation; migrates to a single band at ~87 kDa, observed in Native-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\_002180 (IL-15); NP\_002180.1 (sIL-15R $\alpha$ )

**Tested Applications:** ELISA, Western Blot, CTLL-2 Proliferation Assay

*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 ( $\leq$  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.pblassaysci.com](http://www.pblassaysci.com).

### Authorization

Released by: \_\_\_\_\_ 

Date: September 17, 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.