



Certificate of Analysis

Human IL-15 / sIL-15R α heterodimer, human cell-expressed (HCE)

Catalog No: 11702-1

Lot No:

Size: 10 μ g/vial

Description: Recombinant Human IL-15 / soluble IL-15R α heterodimer, human cell-expressed

Volume: 0.1 ml

Concentration: 0.1 mg/ml

Biological Activity (ED₅₀): <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/ μ g

Molecular Weight: 31 kDa predicted (12.8 kDa for IL-15 + 18.4 kDa for sIL-15R α). 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 α)

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.

Authorization

Released by: _____

Date:

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