

# Recombinant mouse IL-12

CHO expressed mouse interleukin-12 with HSA

Catalog # rcyc-mil12

For research use only, not for diagnostic or therapeutic use

Version # 16C02-MM

## PRODUCT INFORMATION

### Content:

- 10 µg of recombinant mouse IL-12 provided as a white lyophilized powder
- 1.5 ml endotoxin-free water

### Storage and stability:

- Recombinant mouse IL-12 is shipped at room temperature. Upon receipt it should be stored at -20 °C.
- Upon resuspension, prepare aliquots of recombinant mouse IL-12 and store at 4 °C for 1 week or at -20 °C to -80 °C for 12 months.

*Note: Avoid repeated freeze-thaw cycles.*

### Quality control:

- Purity greater than 90% as determined by SDS-PAGE
- Endotoxin: Less than 0.1 EU/µg as determined by the LAL method.
- The biological activity has been confirmed using HEK-Blue™ IL-12 cells (see validation data sheet available on our website).

## DESCRIPTION

Interleukin-12 (IL-12), also known as IL-12p70, cytotoxic lymphocyte maturation factor (CLMF) and natural killer cell stimulatory factor (NKSF), is primarily produced by dendritic cells and macrophages in response to bacterial pathogens<sup>1,2</sup>. It promotes Th1 responses and is a powerful inducer of interferon-gamma (IFN-γ). IL-12 is a heterodimeric 70 kDa cytokine composed of two subunits, p35 and p40.

Recombinant mouse IL-12 is intended for use in cell culture applications. It is produced in Chinese hamster ovary (CHO) cells transfected with a fusion gene encoding the p35 and p40 subunits linked by a decapeptide containing two elastin motifs. These transfected cells produce and secrete biologically active IL-12. Recombinant mouse IL-12 is purified by size-exclusion chromatography.

1. **Vignali DA. & Kuchroo VK., 2012.** IL-12 family cytokines: immunological playmakers. *Nat Immunol.* 13(8):722-8. 2. **Trinchieri G., 1995.** Interleukin-12: a proinflammatory cytokine with immunoregulatory functions that bridge innate resistance and antigen-specific adaptive immunity. *Annu Rev Immunol.* 13:251-76.

## CHARACTERISTICS

**Source:** Mammalian; Chinese hamster ovary (CHO) cells

**Gene Name:** IL-12A\_mouse (subunit p35)

**Gene ID:** 16159

**UniProt ID:** P43431

**Gene Name:** IL-12B\_mouse (subunit p40)

**Gene ID:** 16160

**UniProt ID:** P43432

**Molecular mass:** ≈ 70 kDa

**Formulation:** Recombinant mouse IL-12 was lyophilized from a 0.2 µm filtered phosphate buffer solution (pH 7.4) containing human serum albumin (HSA).

**Solubility:** 100 µg/ml in water

## METHOD

### Preparation of stock solution (100 µg/ml):

1. Add 100 µl endotoxin-free water (provided) to 10 µg of recombinant mouse IL-12.
2. Mix by pipetting. Do not vortex.
3. Prepare aliquots of recombinant mouse IL-12 and store at 4 °C for 1 week or at -20 °C to -80 °C for 12 months. Avoid freeze-thaw cycles.
4. Further dilutions can be prepared in the appropriate aqueous buffer, such as cell culture medium containing serum.

## RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™ IL-12 Cells	hkb-il12

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3-622-34-80

E-mail: [info@invivogen.com](mailto:info@invivogen.com)