

# PRODUCT INFORMATION



## IFN- $\alpha$ Receptor Recognition Peptide 1 (trifluoroacetate salt)

Item No. 44045

**Formal Name:** L-cysteinyl-L-leucyl-L-lysyl-L- $\alpha$ -aspartyl-L-arginyl-L-histidyl-L-aspartic acid, trifluoroacetate salt

**Synonyms:** Cys-Leu-Lys-Asp-Arg-His-Asp-OH, IFN- $\alpha$  (52-58) (human), IFNAR Recognition Peptide 1, IRRP1

**Peptide Sequence:** CLKDRHD-OH

**MF:** C<sub>35</sub>H<sub>59</sub>N<sub>13</sub>O<sub>12</sub>S • XCF<sub>3</sub>COOH

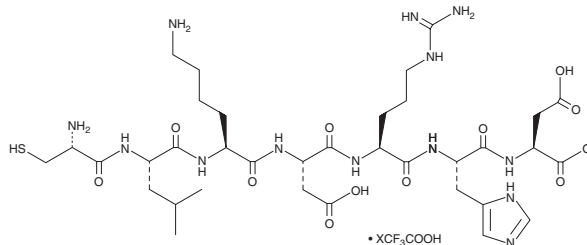
**FW:** 886.0

**Purity:**  $\geq 95\%$

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

IFN- $\alpha$  receptor recognition peptide 1 (IRRP1) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the IRRP1 in water. IRRP1 is slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

IRRP1 is a peptide that corresponds to amino acids 52-58 of full-length IFN- $\alpha$  and is a positive modulator of IFN- $\alpha/\beta$  receptor 1 (IFNAR).<sup>1</sup> It increases the binding capacity of Daudi cells for IFN- $\alpha 2$ . IRRP1 (1  $\mu\text{g}/\text{ml}$ ) enhances the antiproliferative effect of IFN- $\alpha$  in T98G glial cells and enhances the antiviral effect of IFN- $\alpha$  in encephalomyocarditis virus-infected T98G cells.

### Reference

1. Fish, E.N. Interferon receptor recognition peptides enhance the biological potency of interferon alphas. *FEBS Lett.* **365(1)**, 87-91 (1995).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/19/2026

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM