

PRODUCT INFORMATION



KYL (trifluoroacetate salt)

Item No. 36911

Formal Name: L-lysyl-L-tyrosyl-L-leucyl-L-prolyl-L-tyrosyl-L-tryptophyl-L-prolyl-L-valyl-L-leucyl-L-seryl-L-seryl-L-leucine, trifluoroacetate salt

Synonym: KYL Peptide

Peptide Sequence: KYLPYWVPLSSL

MF: $C_{74}H_{108}N_{14}O_{17} \cdot XCF_3COOH$

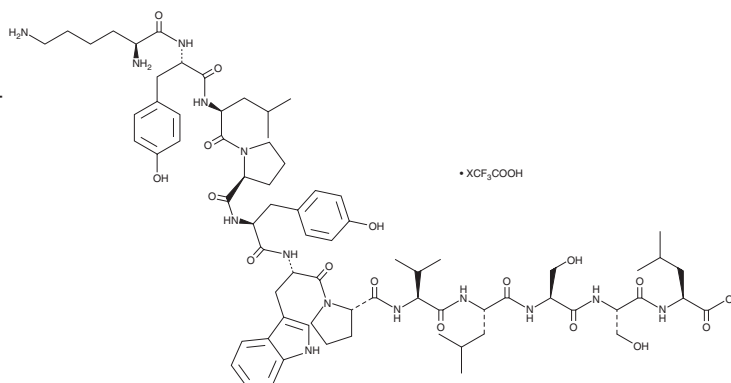
FW: 1,465.7

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

KYL (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the KYL (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. KYL (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KYL (trifluoroacetate salt) in these solvents is approximately 14 and 33 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of KYL (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of KYL (trifluoroacetate salt) in PBS (pH 7.2) is approximately 0.30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

KYL is a 12-amino acid synthetic peptide and an inhibitor of the Eph tyrosine kinase receptor A4 (EphA4; $IC_{50} = 4.22 \mu M$).¹ KYL inhibits EphA4-induced growth cone collapse in embryonic chicken retina explants when used at a concentration of $5 \mu M$.²

References

1. Wu, B., Zhang, Z., Noberini, R., *et al.* HTS by NMR of combinatorial libraries: A fragment-based approach to ligand discovery. *Chem. Biol.* **20**(1), 19-33 (2013).
2. Noberini, R., Koolpe, M., Peddibhotla, S., *et al.* Small molecules can selectively inhibit ephrin binding to the EphA4 and EphA2 receptors. *The Journal of Biological Chemistry* **283**(43), 29461-29472 (2008).

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, 12/21/2022

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