

PRODUCT INFORMATION



GluR2 (831-842) (human, mouse, rat) (trifluoroacetate salt)

Item No. 36966

Formal Name:	L-isoleucyl-L-asparaginy-L-prolyl-L-seryl-L-seryl-L-seryl-L-glutaminy-L-asparaginy-L-seryl-L-glutaminy-L-asparaginy-L-phenylalanine, trifluoroacetate salt	
Synonyms:	AMPA 2, AMPA-selective Glutamate Receptor 2, GluA2 (831-842), Glutamate Ionotropic Receptor AMPA Type Subunit 2, GRIA2 (831-842)	
Peptide Sequence:	INPSSSQNSQNF-OH	H-Ile-Asn-Pro-Ser-Ser-Ser-Gln-Asn-Ser-Gln-Asn-Phe-OH
MF:	$C_{54}H_{83}N_{17}O_{22} \cdot XCF_3COOH$	$\bullet XCF_3COOH$
FW:	1,322.4	
Purity:	≥95%	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

GluR2 (831-842) (human, mouse, rat) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the GluR2 (831-842) (human, mouse, rat) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. GluR2 (831-842) (human, mouse, rat) (trifluoroacetate salt) is slightly soluble in DMSO and dimethyl formamide.

Description

GluR2 (831-842) is a peptide that corresponds to amino acids 831-842 of the C-terminal tail of the ionotropic glutamate receptor subunit GluR2, which is involved in excitatory synaptic transmission.

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, 10/09/2023

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