

PRODUCT INFORMATION



PAR3 (1-6) (human)

Item No. 27129

CAS Registry No.: 320347-28-2

Formal Name: L-threonyl-L-phenylalanyl-L-arginylglycyl-L-alanyl-L-proline

Synonym: TFRGAP

MF: $C_{29}H_{45}N_9O_8$

H—Thr—Phe—Arg—Gly—Ala—Pro—OH

FW: 647.7

Purity: $\geq 98\%$

Supplied as: A crystalline 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

PAR3 (1-6) (human) is supplied as a crystalline solid. A stock solution may be made by dissolving the PAR3 (1-6) (human) in the solvent of choice, which should be purged with an inert gas. PAR3 (1-6) (human) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PAR3 (1-6) (human) in these solvents is approximately 50 mg/ml.

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 PAR3 (1-6) (human) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of PAR3 (1-6) (human) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PAR3 (1-6) is a synthetic peptide agonist of proteinase-activated receptor 1 (PAR1) that corresponds to residues 1-6 of the amino terminal tethered ligand sequence of human PAR3 and residues 39-44 of the full-length human sequence. PAR3 (1-6) activates p42/44 MAPK signaling in fibroblasts expressing PAR1, but not PAR3, an effect that can be blocked by the PAR1 antagonist RWJ 56110.¹

Reference

1. Kaufmann, R., Schulze, B., Krause, G., *et al.* Proteinase-activated receptors (PARs) - The PAR₃ neo-N-terminal peptide TFRGAP interacts with PAR₁. *Regul. Pept.* **125(1-3)**, 61-66 (2005).

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, 11/30/2022

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