# **PRODUCT** INFORMATION



## PBP 10 (trifluoroacetate salt)

Item No. 34535

Formal Name:	N <sup>2</sup> -[2-[3,6- <i>bis</i> (diethylamino) xanthylium-9-yl]benzoyl]-L- glutaminyl-L-arginyl-L-leucyl- L-phenylalanyl-L-glutaminyl-L- valyl-L-lysylglycyl-L-arginine, trifluoroacetate salt	H <sub>2</sub> N NH HN HO I H HO I H HO I H HO H H HO H H H H H H H H H H H H H H
Synonym:	RhB-QRLFQVKGRR-OH	
MF:	C <sub>84</sub> H <sub>127</sub> N <sub>24</sub> O <sub>15</sub> • XCF <sub>3</sub> COOH	
FW:	1,713.1	H <sub>2</sub> N NH NH <sub>2</sub>
Purity:	≥90%	H_N_II
UV/Vis.:	λ <sub>max</sub> : 557 nm	
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

PBP 10 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PBP 10 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. PBP 10 (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 5 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 PBP 10 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of PBP 10 (trifluoroacetate salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

#### Description

PBP 10 is a peptide formyl peptide receptor 2 (FPR2) antagonist (IC<sub>50</sub> = 60 nM).<sup>1</sup> It is active against *E. coli*, *P. aeruginosa*, and *S. pneumoniae* when used at concentrations ranging from 0.1 to 100  $\mu$ M.<sup>2</sup>

#### References

- 1. Skovbakke, S.L., Holdfeldt, A., Nielsen, C., et al. Combining elements from two antagonists of formyl peptide receptor 2 generates more potent peptidomimetic antagonists. J. Med. Chem. 60(16), 6991-6997 (2017).
- 2. Bucki, R., Pastore, J.J., Randhawa, P., et al. Antibacterial activities of rhodamine B-conjugated gelsolin-derived peptides compared to those of the antimicrobial peptides cathelicidin LL37, magainin II, and melittin. Antimicrob. Agents Chemother. 48(5), 1526-1533 (2004).

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

#### SAFFTY 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

uyer 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, 12/05/2022

### 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