

PRODUCT INFORMATION



Thrombin-Sensitive Peptide (ammonium acetate salt)

Item No. 37557

Synonyms:	Click Tag™ Thrombin-Sensitive Peptide, ThS-P
Peptide Sequence:	azidoacetyl alanine-K(5FAM) Azidoacetyl Alanine—Lys—5-FAM—Gly—Ala—Leu—Val—Pro—Arg—Gly—Ser—Ala— GALVPRGSAGK(CPQ2)-NH ₂ Gly—Lys(CPQ2)—NH ₂
FW:	2,143.0
Purity:	≥95% • XC ₂ H ₇ NO ₂
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

Thrombin-sensitive peptide (ThS-P) (ammonium acetate salt) is supplied as a solid. A stock solution may be made by dissolving the ThS-P (ammonium acetate salt) in water. The solubility of ThS-P (ammonium acetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

ThS-P is a synthetic peptide that contains a thrombin-releasable quencher.¹ It has been conjugated to an anti-CD41 antibody *via* click chemistry for use as a platelet-targeting sensor in a microfluidic model of thrombosis and a mouse model of thrombosis to detect increases in fluorescence, which indicate increased thrombin activity.

Reference

1. Welsh, J.D., Colace, T.V., Muthard, R.W., *et al.* Platelet-targeting sensor reveals thrombin gradients within blood clots forming in microfluidic assays and in mouse. *J. Thromb. Haemost.* **10(11)**, 2344-2353 (2012).

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, 08/24/2023

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