

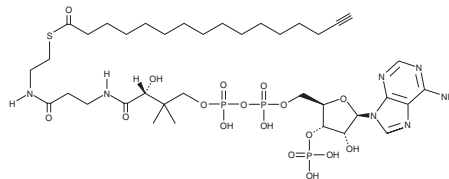
# PRODUCT INFORMATION



## Palmitoyl Alkyne-Coenzyme A

Item No. 36470

**CAS Registry No.:** 1941173-52-9  
**Formal Name:** coenzyme A, S-15-hexadecanoate  
**Synonym:** Click Tag™ Palmitoyl Alkyne-Coenzyme A, Palmitoyl Alkyne-CoA  
**MF:** C<sub>37</sub>H<sub>62</sub>N<sub>7</sub>O<sub>17</sub>P<sub>3</sub>S  
**FW:** 1,001.9  
**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

Palmitoyl alkyne-coenzyme A (palmitoyl alkyne-CoA) is supplied as a solid. A stock solution may be made by dissolving the palmitoyl alkyne-CoA in the solvent of choice, which should be purged with an inert gas. Palmitoyl alkyne-CoA is soluble in organic solvents such as methanol. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Palmitoyl alkyne-CoA is a thioester of  $\omega$ -alkynylated palmitic acid (Item No. 13266) and CoA (Item Nos. 16147 | 21499 | 21722). It has been used as a chemical reporter to label palmitoylated proteins via click chemistry using various probes, including biotin-azide (Item No. 13040) or 3-azido-7-hydroxycoumarin, *in vitro* and *in vivo*.<sup>1</sup>

### Reference

1. Yap, M.C., Kostiuk, M.A., Martin, D.D.O., *et al.* Rapid and selective detection of fatty acylated proteins using  $\omega$ -alkynyl-fatty acids and click chemistry. *J. Lipid Res.* **51**(6), 1566-1580 (2010).

#### 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, 05/08/2024

#### 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](http://WWW.CAYMANCHEM.COM)