

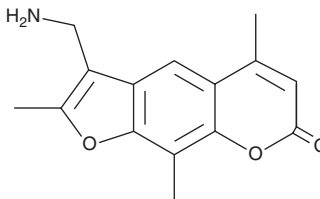
# PRODUCT INFORMATION



## 4'-Aminomethyl-4,5',8-trimethylpsoralen

Item No. 17162

<b>CAS Registry No.:</b>	64358-50-5
<b>Formal Name:</b>	3-(aminomethyl)-2,5,9-trimethyl-7H-furo[3,2-g][1]benzopyran-7-one
<b>Synonym:</b>	4'-Aminomethyltrioxsalen
<b>MF:</b>	C <sub>15</sub> H <sub>15</sub> NO <sub>3</sub>
<b>FW:</b>	257.3
<b>Purity:</b>	≥98%
<b>UV/Vis.:</b>	λ <sub>max</sub> : 240, 287 nm
<b>Supplied as:</b>	A crystalline solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years



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

### Laboratory Procedures

4'-Aminomethyl-4,5',8-trimethylpsoralen is supplied as a crystalline solid. A stock solution may be made by dissolving the 4'-aminomethyl-4,5',8-trimethylpsoralen in the solvent of choice, which should be purged with an inert gas. 4'-Aminomethyl-4,5',8-trimethylpsoralen is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4'-aminomethyl-4,5',8-trimethylpsoralen in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

4'-Aminomethyl-4,5',8-trimethylpsoralen is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4'-aminomethyl-4,5',8-trimethylpsoralen should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 4'-Aminomethyl-4,5',8-trimethylpsoralen has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

4'-Aminomethyl-4,5',8-trimethylpsoralen is a bifunctional photoreactive agent for crosslinking nucleic acids.<sup>1,2</sup> In the dark, it intercalates into double-stranded regions in DNA or RNA. Upon long wavelength UV irradiation (e.g., 365 nm), 4'-aminomethyl-4,5',8-trimethylpsoralen first forms a covalent monoadduct with a pyrimidine base, followed by covalently linking with an adjacent pyrimidine on the opposite strand.<sup>1,2</sup> Short wavelength UV irradiation (e.g., 254 nm) reverses this reaction. 4'-Aminomethyl-4,5',8-trimethylpsoralen has a high binding affinity for RNA and has been used to study different types of RNA from diverse organisms, including viruses.<sup>3-5</sup>

### References

1. Liarakos, C.D., Reinhart, G., Kopper, R.A., *et al.* Characterization of the effects on ovalbumin mRNA of aminomethyl-trimethylpsoralen photoreaction with hen oviduct mRNA. *Nucleic Acids Res.* **15(20)**, 8417-8438 (1987).
2. Garrett-Wheeler, E., Lockard, R.E., and Kumar, A. Mapping of psoralen cross-linked nucleotides in RNA. *Nucleic Acids Res.* **12(7)**, 3405-3423 (1984).
3. Hearst, J.E. and Thiry, L. The photoinactivation of an RNA animal virus, vesicular stomatitis virus, with the aid of newly synthesized psoralen derivatives. *Nucleic Acids Res.* **4(5)**, 1339-1347 (1977).
4. Hui, C.F. and Cantor, C.R. Mapping the location of psoralen crosslinks on RNA by mung bean nuclease sensitivity of RNA-DNA hybrids. *Proc. Natl. Acad. Sci. USA* **82(5)**, 1381-1385 (1985).
5. Raviprakash, K., Sun, P., Raviv, Y., *et al.* Dengue virus photo-inactivated in presence of 1,5-iodonaphthylazide (INA) or AMT, a psoralen compound (4'-aminomethyl-trioxsalen) is highly immunogenic in mice. *Hum. Vaccin. Immunother.* **9(11)**, 2336-2341 (2013).

#### 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, 10/13/2022

#### 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