

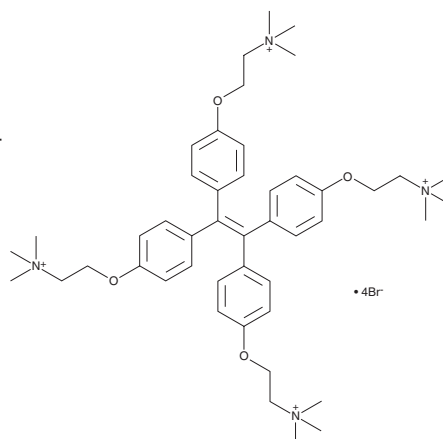
PRODUCT INFORMATION



TTAPE-Me

Item No. 36068

CAS Registry No.: 1219133-97-7
Formal Name: 2,2',2'',2'''-[1,2-ethenediylidenetetrakis(4,1-phenyleneoxy)]tetrakis[N,N,N-trimethylethanaminium, tetrabromide
Synonym: 1,1,2,2-tetrakis[4-(2-Trimethylammonioethoxy)-phenylethene
MF: $C_{46}H_{68}N_4O_4 \bullet 4Br$
FW: 1,060.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 256 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

TTAPE-Me is supplied as a crystalline solid. Aqueous solutions of TTAPE-Me can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of TTAPE-Me in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

TTAPE-Me is an aggregation-induced emission (AIE) fluorogen.¹⁻³ It has been used to detect anionic phospholipids, including cardiolipin, phosphatidylglycerol (PG), and phosphatidylserine (PS), nucleic acids, and the anionic polysaccharide heparin.

References

1. Pyrshev, K., Yesylevskyy, S., and Bogdanov, M. TTAPE-Me dye is not selective to cardiolipin and binds to common anionic phospholipids nonspecifically. *Biophys. J.* **120**(17), 3776-3786 (2021).
2. Hong, Y., Chen, S., Leung, C.W.T., et al. Water-soluble tetraphenylethene derivatives as fluorescent "light-up" probes for nucleic acid detection and their applications in cell imaging. *Chem. Asian J.* **8**(8), 1806-1812 (2013).
3. Wang, Y.J., Lin, L., Zhang, X., et al. Selective, switchable fluorescent probe for heparin based on aggregation-induced emission. *Anal. Biochem.* **514**, 48-54 (2016).

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, 05/30/2023

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