

PRODUCT INFORMATION



LAP

Item No. 43424

CAS Registry No.: 85073-19-4
Formal Name: P-phenyl-P-(2,4,6-trimethylbenzoyl)-
phosphinic acid, monolithium salt

Synonyms: Lithium phenyl-2,4,6-
trimethylbenzoylphosphinate,
Photoinitiator LAP

MF: $C_{16}H_{16}O_3P \cdot Li$

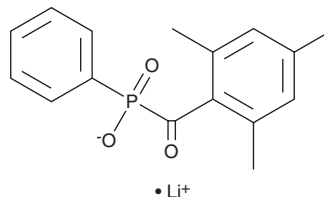
FW: 294.2

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

LAP is supplied as a solid. Aqueous solutions of LAP can be prepared by directly dissolving the solid in aqueous buffers. LAP is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

LAP is a photoinitiator.^{1,2} Upon UV activation, LAP dissociates into benzoyl radicals, which induce polymerization. LAP has been used to induce polymerization and hardening of methacrylated hydrogels for vascular tissue engineering and the development of stem cell spheroids.

References

1. Martins, E.A.G., Deus, I.A., Gomes, M.C., *et al.* Human chorionic membrane-derived tunable hydrogels for vascular tissue engineering strategies. *Adv. Healthc. Mater.* **14(1)**, e2401510 (2025).
2. Zhu, S., Liao, X., Xu, Y., *et al.* 3D bioprinting of high-performance hydrogel with in-situ birth of stem cell spheroids. *Bioact. Mater.* **43**, 392-405 (2024).

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, 06/17/2025

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