

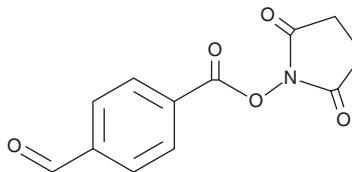
# PRODUCT INFORMATION



## Succinimidyl 4-formylbenzoate

Item No. 30477

**CAS Registry No.:** 60444-78-2  
**Formal Name:** 4-formyl-benzoic acid, 2,5-dioxo-1-pyrrolidinyl ester  
**Synonyms:** SFB, Succinimidyl *p*-formylbenzoate  
**MF:** C<sub>12</sub>H<sub>9</sub>NO<sub>5</sub>  
**FW:** 247.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 253 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

Succinimidyl 4-formylbenzoate is supplied as a crystalline solid. A stock solution may be made by dissolving the succinimidyl 4-formylbenzoate in the solvent of choice, which should be purged with an inert gas. Succinimidyl 4-formylbenzoate is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of succinimidyl 4-formylbenzoate in these solvents is approximately 30 mg/ml.

Succinimidyl 4-formylbenzoate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, succinimidyl 4-formylbenzoate should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Succinimidyl 4-formylbenzoate has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Succinimidyl 4-formylbenzoate is a heterobifunctional building block.<sup>1,2</sup> It has been used in the conjugation of oligonucleotides to antibodies for protein detection, as well as in the synthesis of fluorescent BODIPY-etoposide.

### References

1. Kozlov, I.A., Melnyk, P.C., Stromborg, K.E., *et al.* Efficient strategies for the conjugation of oligonucleotides to antibodies enabling highly sensitive protein detection. *Biopolymers* **73(5)**, 621-630 (2004).
2. Lazo, J.S., Li, T., Woo, E.S., *et al.* Chemical synthesis and biological activity of a novel fluorescent etoposide derivative. *Biochem. Pharmacol.* **53(5)**, 15-22 (1997).

#### 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, 05/13/2020

#### 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