

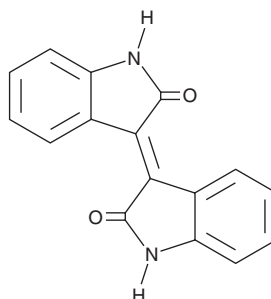
PRODUCT INFORMATION



Isoindigo

Item No. 33208

CAS Registry No.: 476-34-6
Formal Name: 3-(1,2-dihydro-2-oxo-3H-indol-3-ylidene)-1,3-dihydro-2H-indol-2-one
Synonym: Isoindigotin
MF: C₁₆H₁₀N₂O₂
FW: 262.3
Purity: ≥98%
UV/Vis.: λ_{max}: 267 nm
Supplied as: A crystalline 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

Isoindigo is supplied as a crystalline solid. A stock solution may be made by dissolving the isoindigo in the solvent of choice, which should be purged with an inert gas. Isoindigo is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of isoindigo in these solvents is approximately 10 mg/ml.

Description

Isoindigo is an electron-accepting building block.^{1,2} It has been used in the synthesis of polymers for use in organic photovoltaics.

References

1. Lei, T., Wang, J.-W. and Pei, J. Design, synthesis, and structure-property relationships of isoindigo-based conjugated polymers. *Acc. Chem. Res.* **47**(4), 1117-1126 (2014).
2. Stalder, R., Mei, J., Graham, K.R., *et al.* Isoindigo, a versatile electron-deficient unit for high-performance organic electronics. *Chem. Mater.* **26**(1), 664-678 (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, 11/03/2022

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