

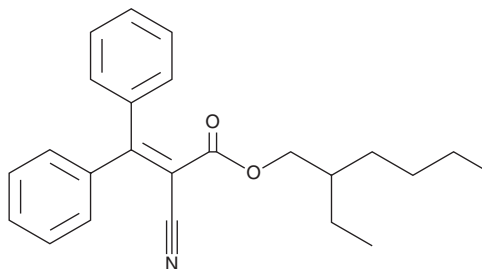
PRODUCT INFORMATION



Octocrylene

Item No. 35461

CAS Registry No.: 6197-30-4
Formal Name: 2-cyano-3,3-diphenyl-2-propenoic acid
2-ethylhexyl ester
Synonym: 2-Ethylhexyl 2-cyano-3,3-diphenylacrylate
MF: $C_{24}H_{27}NO_2$
FW: 361.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 303 nm
Supplied as: A liquid
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

Octocrylene is supplied as a liquid. A stock solution may be made by dissolving the octocrylene in the solvent of choice, which should be purged with an inert gas. Octocrylene is soluble in the organic solvent DMSO.

Description

Octocrylene is a UV filter.^{1,2} It stabilizes the full-spectrum UVA blocker avobenzone (Item No. 23836) to UV radiation.¹ Octocrylene inhibits 17β -estradiol- or dihydrotestosterone-induced transactivation of estrogen receptor α (ER α) or the androgen receptor in *S. cerevisiae* expressing the human receptors (IC_{50} s = 2.17 and 0.627 mM, respectively).² Topical application of octocrylene (15 or 30% w/v) increases thymidine incorporation in a mouse local lymph node assay (LLNA), indicating allergenic activity.³ Octocrylene has been found in influent and effluent wastewater.⁴ Formulations containing octocrylene have been used in sunscreen products.

References

1. Beasley, D.G. and Meyer, T.A. Characterization of the UVA protection provided by avobenzone, zinc oxide, and titanium dioxide in broad-spectrum sunscreen products. *Am. J. Clin. Dermatol.* **11**(6), 413-421 (2010).
2. Kunz, P.Y. and Fent, K. Multiple hormonal activities of UV filters and comparison of *in vivo* and *in vitro* estrogenic activity of ethyl-4-aminobenzoate in fish. *Aquat. Toxicol.* **79**(4), 305-324 (2006).
3. Karlsson, I., Broecke, K.V., Mårtensson, J., et al. Clinical and experimental studies of octocrylene's allergenic potency. *Contact Dermatitis* **64**(6), 343-352 (2011).
4. Balmer, M., Buser, H.-R., Müller, M.D., et al. Occurrence of some organic UV filters in wastewater, in surface waters, and in fish from Swiss lakes. *Environ. Sci. Technol.* **39**(4), 953-962 (2005).

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, 02/18/2026

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