

PRODUCT INFORMATION

N-[(3-(Anilinomethylene)-2-chloro-1-cyclohexen-1-yl)methylene]aniline (hydrochloride)

Item No. 30560

CAS Registry No.: 63857-00-1

Formal Name: N-[[2-chloro-3-[(phenylamino)methylene]-1-cyclohexen-1-yl]methylene]benzenamine, monohydrochloride

MF: C₂₀H₁₉ClN₂ • HCl

FW: 359.3

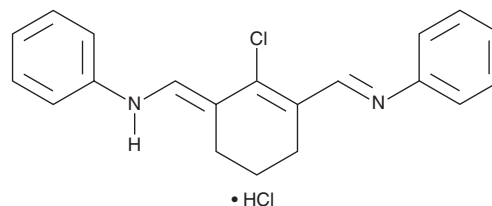
Purity: ≥98%

UV/Vis.: λ_{max}: 523 nm

Supplied as: A 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

N-[(3-(Anilinomethylene)-2-chloro-1-cyclohexen-1-yl)methylene]aniline (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the N-[(3-(anilinomethylene)-2-chloro-1-cyclohexen-1-yl)methylene]aniline (hydrochloride) in the solvent of choice, which should be purged with an inert gas. N-[(3-(Anilinomethylene)-2-chloro-1-cyclohexen-1-yl)methylene]aniline (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N-[(3-(anilinomethylene)-2-chloro-1-cyclohexen-1-yl)methylene]aniline (hydrochloride) in these solvents is approximately 1, 3, and 0.25 mg/ml, respectively.

Description

N-[(3-(Anilinomethylene)-2-chloro-1-cyclohexen-1-yl)methylene]aniline is a building block.^{1,2} It has been used in the synthesis of various cyanine-containing dyes.

References

1. Rivera, L., Puyol, M., Miltsov, S., *et al.* New hexamethine-hemicyanine dyes for the development of integrated optochemical sensors. *Anal. Bioanal. Chem.* **387**(6), 2111-2119 (2007).
2. Henary, M., Pannu, V., Owens, E.A., *et al.* Near infrared active heptacyanine dyes with unique cancer-imaging and cytotoxic properties. *Bioorg. Med. Chem. Lett.* **22**(2), 1242-1246 (2012).

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