

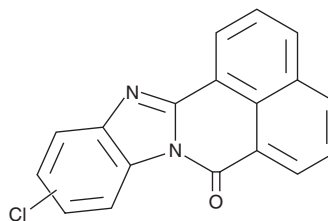
PRODUCT INFORMATION



10(11)-Cl-BBQ

Item No. 29978

Formal Name: 10(or 11)-chloro-7H-benzimidazo[2,1-a]benz[de]isoquinolin-7-one
MF: C₁₈H₉ClN₂O
FW: 304.7
Purity: ≥95% (mixture of isomers)
UV/Vis.: λ_{max}: 219, 243, 285, 296, 382 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

10(11)-Cl-BBQ is supplied as a crystalline solid. A stock solution may be made by dissolving the 10(11)-Cl-BBQ in the solvent of choice, which should be purged with an inert gas. 10(11)-Cl-BBQ is soluble in the organic solvent DMSO.

Description

10(11)-Cl-BBQ contains the 10-Cl-BBQ and 11-Cl-BBQ congeners of Cl-BBQ and acts as an aryl hydrocarbon receptor (AhR) agonist.¹ It induces transcription of the AhR target cytochrome P450 (CYP) isoform *Cyp1a1* in mouse liver when administered intraperitoneally at a dose of 10 mg/kg. This product is a mixture of 10-Cl-BBQ and 11-Cl-BBQ.

Reference

1. Ehrlich, A.K., Pennington, J.M., Bisson, W.H., et al. TCDD, FICZ, and other high affinity AhR ligands dose-dependently determine the fate of CD4⁺ T cell differentiation. *Toxicol. Sci.* **161**(2), 310-320 (2018).

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, 05/02/2023

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