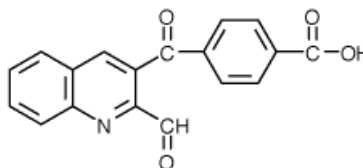


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: CBQCA

CATALOG # 90070

MOLECULAR INFORMATION: C₁₈H₁₁NO₄
Mwt: 305
[131124-59-9]

**PROPERTIES:**

Color & Form Light yellow solid
Solubility Soluble in DMF or DMSO
Absorption/Emission 465/560nm (MeOH) for reaction product with amines

STORAGE AND HANDLING: Store at -20°C and protect from light

APPLICATION: CBQCA (3-(4-carboxybenzoyl)quinoline-2-carboxaldehyde) reacts specifically with primary amines to form conjugates that can be analyzed by electrophoretic or chromatographic methods. CBQCA also reacts with hydrophilic peptides and amino sugars. The amine derivatized product (7-aza-1-cyano-5, 6-benzisindole) is highly fluorescent that is optimally excited at 450 nm and have emission maximum at 560 nm.

Ref.: 1) *Anal. Chem.* **63**, 408 (1991); 2) *Anal. Chem.* **63**, 413 (1991); 3) *J. Chromatography*, **559**, 223 (1991).

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--