

CBQCA

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.



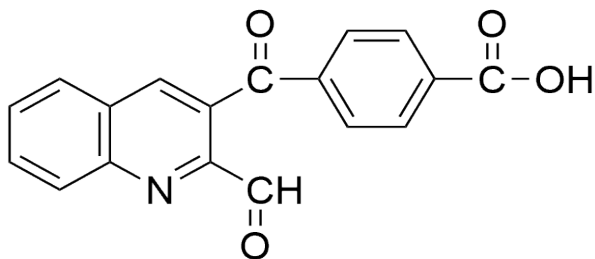
Product attributes

Excitation/Emission	465/560 nm (reaction product)
---------------------	-------------------------------

Product Description

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.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 465/560 \text{ nm}$ (for reaction product with amines)
- Light yellow solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- $\text{C}_{18}\text{H}_{11}\text{NO}_4$
- MW: 305



This datasheet was generated on June 10, 2026 at 10:05:17 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/cbqca/>