## Product attributes

Call us: 800-304-5357

Excitation/Emission

465/560 nm (reaction product)

## **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 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-benzisoindole) is highly fluorescent.

- $\lambda_{Ex}/\lambda_{Em}(MeOH) = 465/560$  nm (for reaction product with amines)
- Light yellow solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- C<sub>18</sub>H<sub>11</sub>NO<sub>4</sub>
- MW: 305

This datasheet was generated on November 30, 2025 at 10:56:32 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cbqca/">https://biotium.com/product/cbqca/</a>