

CBQCA Reagent [3-(4-carboxybenzoyl)quinoline-2-carboxaldehyde]

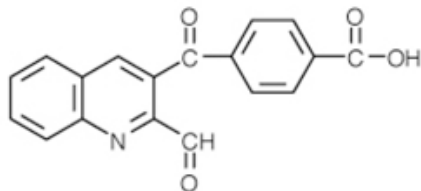
Catalog Number	Packaging Size
C300	25 mg

Storage upon receipt: -20°C, protect from light

Introduction

The **CBQCA reagent** (Known as ATTO-TAG™, Trade Mark of Molecular Probes) reacts specifically with primary amines to form conjugates that can be analyzed by electrophoretic or chromatographic methods. The resulting products of CBQCA are maximally excited at 450 nm and have emission maxima at ~550 nm. In capillary zone electrophoresis (CZE), the sensitivity of detection of CBQCA conjugates should be in the attomole range (10⁻¹⁸ moles). The high sensitivity, freedom from background and long-wavelength excitability make these potential reagents for researcher, diagnostic and forensic applications.

Specifications

Label:	ATTO-TAG	
Ex/Em:	450/550 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₁₈ H ₁₁ NO ₄	
Molecular Weight:	305.29	
CAS Number:	-	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [Mycothiol biosynthesis and metabolism. Cellular levels of potential intermediates in the biosynthesis and degradation of mycothiol in mycobacterium smegmatis.](#)
Anderberg SJ, Newton GL, Fahey RC
J Biol Chem (1998) 273:30391-30397
2. [Capillary electrophoresis laser-induced fluorescence for screening combinatorial peptide libraries in assays of botulinum neurotoxin A.](#)
Laing TD, Marenco AJ, Moore DM, Moore GJ, Mah DC, Lee WE,
J Chromatogr B Analyt Technol Biomed Life Sci (2006) 843:240-246