

MEQ

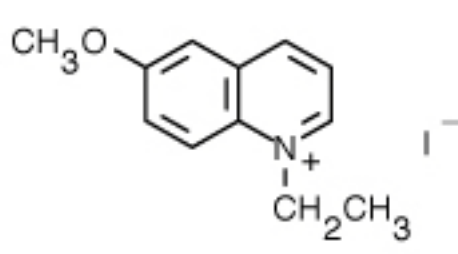
Catalog Number	Packaging Size
C273	50 mg

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

Introduction

The chloride indicator **MEQ** has absorption/emission maxima ~344/440. MEQ can be reduced to create the cell-permeant dihydro-MEQ (DiH-MEQ) which can be passively loaded into cells and subsequently oxidized to the chloride-sensitive MEQ.

Specifications

Label:	MEQ	
Ex/Em:	344/440 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula	C ₁₂ H ₁₄ INO	
Molecular Weight:	315.15	
CAS Number:	-	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Chloride indicator

References:

Glycosylphosphatidylinositol-anchored proteins and actin cytoskeleton modulate chloride transport by channels formed by the *Helicobacter pylori* vacuolating cytotoxin VacA in HeLa cells.

Gauthier NC, Ricci V, Gounon P, Doye A, Tauc M, Poujeol P, Boquet P
J Biol Chem (2004) 279:9481-9489

Cell swelling stimulates cytosol to membrane transposition of ICln.

Ritter M, Ravasio A, Jakab M, Chwatal S, Fürst J, Laich A, Gschwentner M, Signorelli S, Burtscher C, Eichmüller S, Paulmichl M
J Biol Chem (2003) 278:50163-50174