

# Chloride Channel Assay Kit (Colorimetric) ab176767

★★★★★ [1 Abreviews](#) [2 References](#) [1 Image](#)

## Overview

Product name	Chloride Channel Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Adherent cells, Suspension cells
Assay type	Quantitative
Product overview	<p>Abcam's Chloride Channel Assay Kit (Colorimetric) (ab176767) provides a sensitive and robust colorimetric method for studying chloride channels. The assay is based on our proprietary iodide indicator (Iodide Sensor Blue dye) to measure iodide concentration, as low as 30 nM of iodide can be detected. Iodide Sensor Blue dye forms a blue complex with iodide, which has absorption spanning from the UV to 700 nm. Thus a few absorption wavelengths can be used for monitoring the iodide-dependent color change.</p> <p>Abcam's Chloride Channel Assay Kit (Colorimetric) (ab176767) provides an optimized assay method for monitoring chloride channels. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format.</p>

Notes	<p>Chloride channels have a variety of important physiological and cellular functions that include regulation of pH, volume homeostasis, organic solute transport, cell migration, cell proliferation and differentiation. Chloride channels represent valuable drug targets. A number of chronic disease states such as cystic fibrosis and Bartter's syndrome are due to defects in chloride channel functions. However, the existing technologies for screening chloride channel modulators are a compromise between throughput, sensitivity and physiological relevance.</p> <p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p>
-------	--

Platform	Microplate reader
----------	-------------------

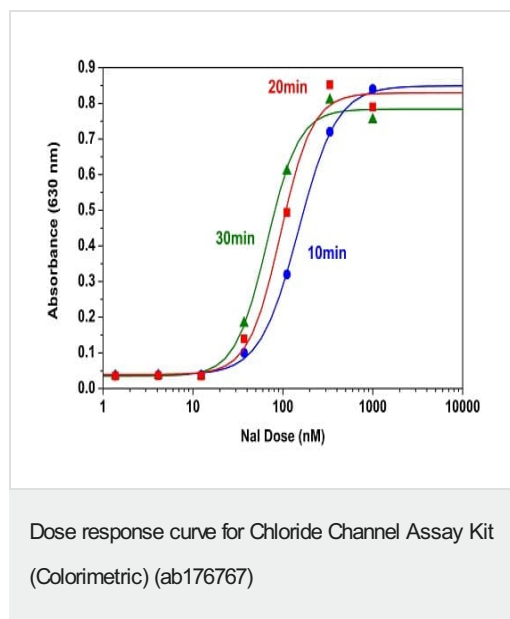
## Properties

Storage instructions	Store at +4°C. Please refer to protocols.
----------------------	---

Components	10 x 96 tests
100X Iodide Sensor Enhancer	1 x 0.5ml

Components	10 x 96 tests
10X Cell Lysis Buffer	1 x 5ml
I-Loading Buffer	1 x 100ml
Iodide Sensor Blue dye	1 x 50ml

## Images



Sodium iodide (NaI) dose response was measured with Abcam's Chloride Channel Assay Kit (Colorimetric) (ab176767) in a 96-well black wall/clear bottom plate. As low as 30 nM of NaI was detected with 10 minutes incubation time (n = 3).

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors