

Amplite™ Colorimetric Hypochlorite (Hypochlorous Acid) Assay Kit

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 13845 (200 tests)	Frozen	Absorbance microplate reader

Introduction

Hypochlorite anion (ClO^-) and its protonated form, hypochlorous acid (HClO) are critical reactive oxygen species (ROS) in biological systems. Uncontrolled production of hypochlorite (hypochlorous acid) can lead to tissue damage and diseases including arthritis, renal failure and cancers. In addition, sodium hypochlorite (NaClO) has been widely used as a bleaching agent for surface cleaning, odor removal and water disinfection in our daily life. Exposure to large amount of sodium hypochlorite can lead to poisoning with the symptoms of serious breathing problems, stomach irritation, redness and pain on skin and eye. Therefore, highly selective and sensitive detection of hypochlorite (hypochlorous acid) is of toxicological and environmental importance. Amplite™ Colorimetric Hypochlorite (Hypochlorous Acid) Assay Kit offers a sensitive assay for measuring hypochlorite (hypochlorous acid) with high specificity. The Oxirite™ Hypochlorite Sensor selectively reacts with hypochlorite (hypochlorous) to generate a red color product. The assay can be measured with an absorbance microplate reader around 550 nm.

Kit Components

Components	Amount
Component A: Oxirite™ Hypochlorite Sensor (light sensitive)	1 vial
Component B: Assay Buffer	1 bottle (20 mL)
Component C: Hypochlorite Standard	1 vial (300 μL)
Component D: DMSO	1 vial (600 μL)

Assay Protocol for One 96-Well Plate

Brief Summary

Prepare hypochlorite assay mixture (50 μL) → Add hypochlorite or test samples (50 μL) → Incubate at room temperature for 3-5 min → Monitor absorbance increase at OD of 555±5 nm

Note: Thaw one vial of each kit component at room temperature before starting the experiment.

1. Prepare Oxirite™ hypochlorite sensor stock solution (20X):

Add 500 μL of DMSO into the vial of Oxirite™ hypochlorite sensor (**Component A**) to make 20X stock solution.
Note: Make single use aliquots, and store unused 20X Oxirite™ hypochlorite sensor stock solution at -20°C , avoid light and repeat freeze-thaw cycles.

2. Prepare hypochlorite assay mixture:

Add 250 μL of Oxirite™ hypochlorite sensor stock solution (from Step 1) into 5 mL of assay buffer (**Component B**), and mix well to make hypochlorite assay mixture (**Component A+B**).
Note: This hypochlorite assay mixture is enough for one 96-well plate. It is not stable, use it promptly.

3. Prepare serial dilutions of hypochlorite standards (0 to 1%):

3.1 Prepare hypochlorite standard dilutions: Add 100 μL of hypochlorite standard (**Component C**) into 400 μL of assay buffer (**Component B**) to get 1% hypochlorite standard solution. Perform 1:3 serial dilutions in assay buffer (**Component B**) get approximately 1%, 0.3%, 0.1%, 0.03%, 0.01%, 0.003%, 0.001% and 0% serially diluted hypochlorite standards.

Note: The unused hypochlorite standard (Component C) should be divided into single use aliquots and stored at -20°C .

3.2 Add serial dilutions of hypochlorite standard and hypochlorite/hypochlorous acid containing test samples into a 96-well clear bottom microplate microplate as described in Tables 1 and 2.

