

COX Fluorescent Inhibitor Screening Assay Kit Short Protocol

Item No. 700100

REAGENT PREPARATION

1. **Assay Buffer (10X) (Item No. 760114)** - Dilute 3 mls with 27 ml HPLC-grade water (final formulation is 100 mM Tris-HCl, pH 8.0); stable 6 months at 4°C.
2. **Heme (Item No. 760116)** - Dilute 40 µl with 960 µl diluted assay buffer; stable 12 hours at 22°C.
3. **COX-1 (ovine) (Item No. 700103)** - For approximately 48 wells, dilute 40 µl of enzyme with 440 µl of diluted Assay Buffer and store on ice; use within 1 hour.
4. **COX-2 (human recombinant) (Item No. 700104)** - For approximately 48 wells, dilute 40 µl of enzyme with 440 µl of diluted Assay Buffer and store on ice; use within 1 hour.
5. **Arachidonic Acid (Item No. 760113)** - Transfer 100 µl of the supplied solution to another vial, add 100 µl of KOH (Item No. 760115), vortex, and dilute with 800 µl of HPLC-grade water (concentration = 2 mM); use within 30 minutes.
6. **Potassium Hydroxide (Item No. 760115)** - Ready to use as supplied.
7. **DMSO (Item No. 700001)** - Ready to use as supplied.
8. **ADHP (Item No. 700002)** - Dissolve the contents of one vial with 100 µl DMSO (Item No. 700001) and then add 900 µl of diluted Assay Buffer; use within 30 minutes.
9. **DuP-697 (Item No. 760158)**; optional - Ready to use as supplied. Assaying 10 µl will yield approximately 90% inhibition or greater.
10. **SC-560 (Item No. 760159)**; optional - Ready to use as supplied. Assaying 10 µl will yield approximately 90% inhibition or greater.



Short Protocol
Item 700100

Tech Support
888.526.5351
techserv@caymanchem.com

COX Fluorescent Inhibitor Screening Assay Kit Short Protocol

Item No. 700100

PERFORMING THE ASSAY

Pipette the reagents as indicated in the table below and incubate the reaction as described.

Wells	Assay Buffer	Heme	Enzyme	Compound Solvent	Inhibitor/ Sample
100% Initial Activity (IA) 3 wells	150 µl	10 µl	10 µl	10 µl	-
Background (IA) 3 wells	160 µl	10 µl	-	10 µl	-
Inhibitor/Sample triplicate recommended	150 µl	10 µl	10 µl	-	10 µl
Incubate inhibitor + enzyme for 5 minutes at room temperature					
Add 10 µl ADHP to each of the wells					
Quickly add 10 µl Arachidonic Acid to all wells to initiate reactions					
Incubate precisely 2 minutes at room temperature					
Read the plate at Excitation = 530-540; Emission = 585-595					

CALCULATIONS

$$\% \text{ Inhibition} = \frac{[\text{Initial Activity} - \text{Sample Activity}]}{\text{Initial Activity}} \times 100$$



Short Protocol
Item 700100

Tech Support
888.526.5351
techserv@caymanchem.com

For Research Use Only
www.caymanchem.com

09/02/2015