

PRODUCT: DiscoveryPak™ Stem Cell Fate Regulator Set X**CATALOG #:** K396-2**AMOUNT:** 2 Modulators**STORAGE:** Store at -20°C. Protect from light.**DESCRIPTION:** A convenient set of two small molecules useful for increasing the cloning efficiency in individualized human pluripotent stem cells (hPSCs). The two products are shown in the following table:

Product	Cat. No.	Biological Activity	Size
Pioglitazone	1877-5	A PPAR γ agonist	5 mg
Y-27632, DiHCl	1596-5	A ROCK inhibitor	5 mg

REFERENCE: Kajabadi, N.S., *et al.* (2015). *J. Biol. Chem.* **290**, 26303-13**HANDLING:** Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.**RELATED PRODUCTS:**

DiscoveryPak™ Stem Cell Fate Regulator Set I (**Cat. No. K852-3**)
 DiscoveryPak™ Stem Cell Fate Regulator Set II (**Cat. No. K853-8**)
 DiscoveryPak™ Stem Cell Fate Regulator Set III (**Cat. No. K854-3**)
 DiscoveryPak™ Stem Cell Fate Regulator Set IV (**Cat. No. K855-3**)
 DiscoveryPak™ Stem Cell Fate Regulator Set V (**Cat. No. K865-5**)
 DiscoveryPak™ Stem Cell Fate Regulator Set VI (**Cat. No. K866-5**)
 (±) –Bay K 8644 (**Cat. No. 1682-5**)
 BIX01294 (**Cat. No. 1678-5, 25**)
 CHIR99021 (**Cat. No. 1677-5, 25**)
 EZSolution™ CHIR99021 (**Cat. No. 1748-5**)
 PD0325901 (**Cat. No. 1643-2**)
 PS-48 (**Cat. No. 1869-5**)
 Pyrintegrin (**Cat. No. 1729-1,5**)
 EZSolution™ Pyrintegrin (**Cat. No. 1737-1**)
 RepSox (**Cat. No. 1894-5, 25**)
 Reversine (**Cat. No. 1851-1, 5**)
 RG 108 (**Cat. No. 1679-10,30**)
 SB-431542 (**Cat. No. 1674-1**)
 EZSolution™ SB-431542 (**Cat. No. 1872-1**)
 Sodium Butyrate (**Cat. No. 1609-100, 1000**)
 Thiazovivin (**Cat. No. 1681-1,5**)
 EZSolution™ Thiazovivin (**Cat. No. 1736-1**)
 Tranylcypromine Hemisulfate (Parnate) (**Cat. No. 1816-25,100**)
 Valproic Acid, Sodium Salt (**Cat. No. 1647-200**)
 Y-27632,2 HCl (**Cat. No. 1596-1, 5, 50**)
 EZSolution™ Y-27632, HCl (**Cat. No. 1784-5**)

USAGE: *FOR RESEARCH USE ONLY! Not to be used in humans.*