



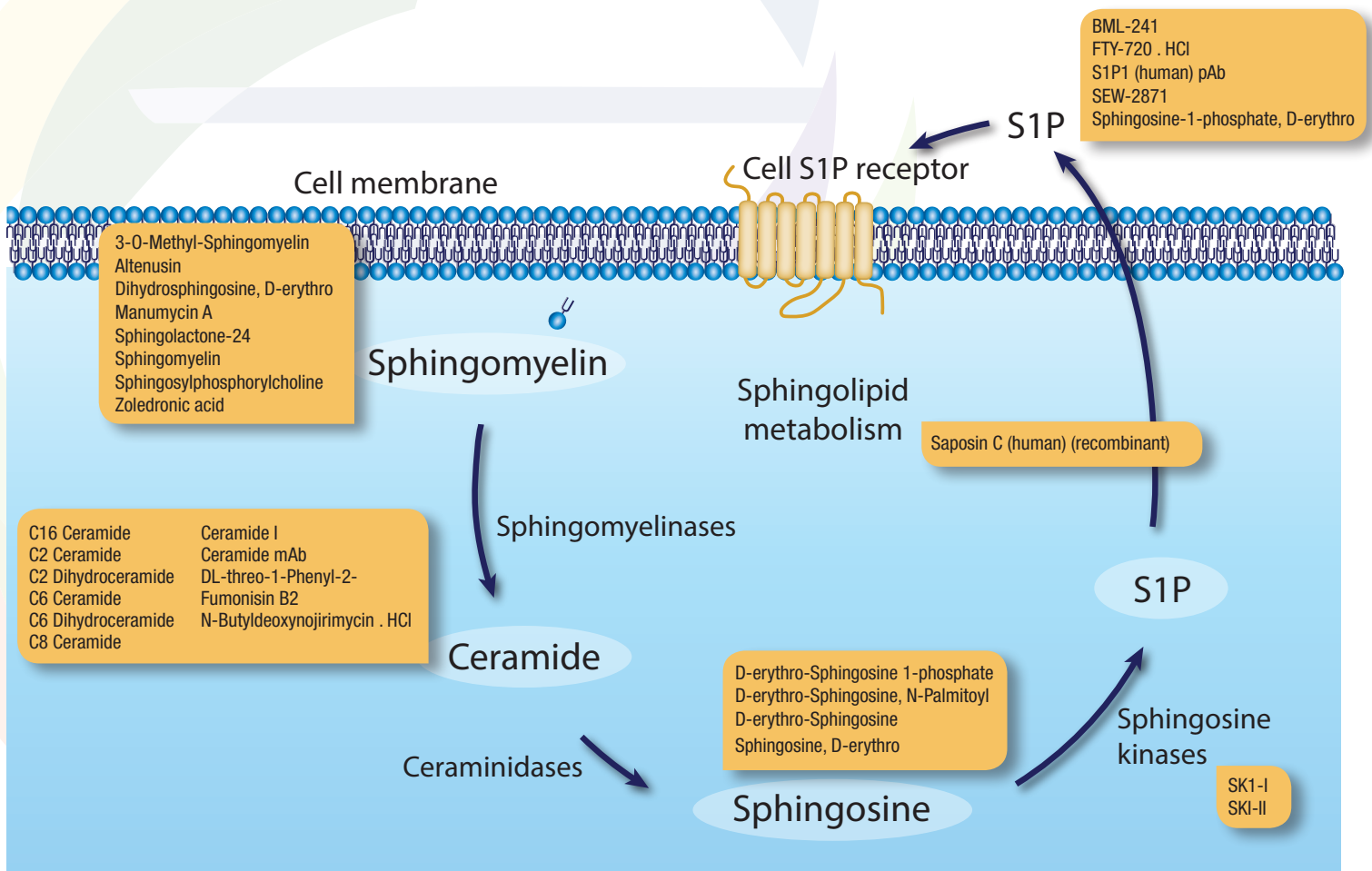
Exclusive Sphingosine Kinase 1 Inhibitor

Sphingosine Kinase 1 Inhibitor (SK1-I)

A sphingosine analog and competitive inhibitor specific for sphingosine kinase 1

Sphingosine Kinase 1 Inhibitor (BML-EI411)

Sphingolipids are a versatile class of membrane lipid. Sphingolipid metabolites, such as ceramide, sphingosine, and sphingosine-1-phosphate are lipid signaling molecules which are involved in diverse cellular processes. Sphingosine can be phosphorylated *in vivo* via two kinases, sphingosine kinase type 1 and sphingosine kinase type 2, leading to the formation of sphingosine-1-phosphate (S1P). Sphingosine-1-phosphate is a potent signaling lipid involved in numerous biological functions including, cellular proliferation, survival, migration, invasion, differentiation, architecture, immune cell trafficking, angiogenesis and vascular integrity.



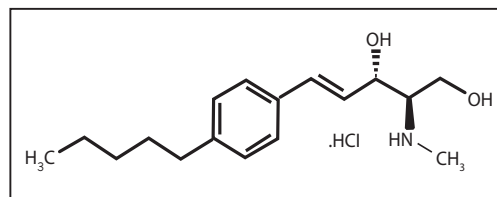
Fuel Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS
 Enzo Life Sciences Inc.
 10 Executive Boulevard
 Farmingdale, NY 11735
 Phone: 800.942.0430
 Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE
 Enzo Life Sciences (ELS) AG
 Industriestrasse 17
 CH-4415 Lausen, Switzerland
 Phone: +41 61 926 8989
 Fax: +41 61 926 8979
info-eu@enzolifesciences.com

RELATED PRODUCTS			
Product Category	Product Name	Product #	
	alpha-Mannosyl Ceramide	ALX-306-031	
Activator	C16 Ceramide	BML-SL115	
	C2 Ceramide	BML-SL100	
	C2-Phytoceramide	BML-SL151	
	C6 Ceramide	BML-SL110	
	C8 Ceramide	BML-SL112	
	C8 ceramine, D-erythro	BML-SL145	
	D-erythro-Sphingosine 1-phosphate (caged)	ALX-306-017	
	D-erythro-Sphingosine, N-Palmitoyl-	ALX-306-006	
	Sphingosine-1-phosphate, D-erythro	BML-SL140	
	Sphingosylphosphorylcholine	BML-SL141	
Antibody	Ceramide , mAb (MID 15B4)	ALX-804-196	
	S1P1 (human), pAb	BML-SA557	
Inactive Control	C2 Dihydroceramide	BML-SL101	
	C6 Dihydroceramide	BML-SL111	
	C8-Dihydroceramide	BML-SL113	
Inhibitor	SK1-I	BML-EI411	
	3-O-Methyl-Sphingomyelin	BML-SL225	
	Altenusin	ALX-350-325	
	D-erythro-Sphingosine (synthetic)	ALX-306-011	
	D-erythro-Sphingosine chloride, N,N,N-Trimethyl-	ALX-306-022	
	Dihydrosphingosine, D-erythro	BML-SL125	
	DL-threo-1-Phenyl-2-decanoylamino-3-morpholino-1-propanol	BML-SL210	
	DL-threo-1-Phenyl-2-palmitoylamino-3-morpholino-1-propanol	BML-SL215	
	Epoxyquinone G109 (racemic)	ALX-270-341	
	Fumonisin B2	BML-SL219	
	Manumycin A	ALX-350-241	
	N-Butyldeoxyojirimycin . HCl	BML-SL230	
	SKI-II	BML-EI406	
	Sphingolactone-24	ALX-270-453	
	Sphingosine, D-erythro	BML-EI155	
	Zoledronic acid	ALX-430-153	
	Ligand	BML-241	BML-SL227
		Dihydrosphingosine-1-phosphate, D-erythro	BML-SL143
		FTY-720 . HCl	BML-SL233
SEW-2871		BML-LP106	
Protein	S1P receptor 1 (rat), (membrane preparation)	BML-SE327	
	Saposin C (human), (recombinant)	ALX-201-262	
	Sphingomyelinase (Staphylococcus aureus)	BML-SE108	
	Sphingosine kinase 1 (human), (recombinant) (His-tag)	BML-SE424	
	Sphingosine kinase 2 (human), (recombinant) (His-tag)	BML-SE425	
Sphingolipid	3-Ketosphingosine	BML-SL130	
	C8 Ceramide-1-phosphate, D-erythro	BML-SL142	
	Ceramide I (bovine brain)	ALX-303-001	
	Ceramide II (bovine brain)	ALX-303-002	
	D-erythro-Sphingosine, N-Stearoyl-	ALX-306-007	
	D-erythro-Sphingosine, N-Stearoyl- (synthetic)	ALX-306-023	
	N,N-Dimethylsphingosine, D-erythro	BML-SL105	
	Phytosphingosine	BML-SL150	
Sphingomyelin	BML-SL135		

Highly Specific Inhibitor for SK1



Sphingosine kinase 1 inhibitor (BML-EI411)

Benefits

- Competitive inhibitor specific for sphingosine kinase 1 (SK1)
- Exclusive to Enzo Life Sciences
- High purity $\geq 98\%$ (HPLC)
- Highly specific (does not inhibit SK2, PKCa, PKCd, PKA, AKT1, ERK1, EGFR, CDK2, IKKb or CamK2b)
- Decreases sphingosine-1-phosphate levels and increases proapoptotic precursor ceramide
- SK1-I inhibits tumor cell growth *in vitro* and *in vivo*

Research Areas

- Atherosclerosis
- Cancer
- Diabetes
- Hepatitis C
- Inflammatory disorders
- Respiratory distress

Citations:

A novel E2F/Sphingosine kinase 1 axis regulates anthracycline response in squamous cell carcinoma: M. Hazar-Rethinam, et al.; Clin. Cancer Res. 21, 417 (2015), **Application(s):** Inhibition of Sphingosine kinase 1 in doxorubicin-treated SCC cells and *in vivo*.

Inhibition of Sphingosine Kinase 1 Ameliorates Angiotensin II-induced Hypertension and Inhibits Transmembrane Calcium Entry via Store-Operated Calcium Channel: P. C. Wilson, et al.; Mol. Endocrinol. (2015), **Application(s):** Cell Culture.

Sphingosine Kinases Signalling in Carcinogenesis: G. Marfe, et al.; Mini Rev. Med. Chem. 15, 300 (2015), **Application(s):** Inhibition of Sphingosine kinase 1.

CONTACT INFORMATION

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com