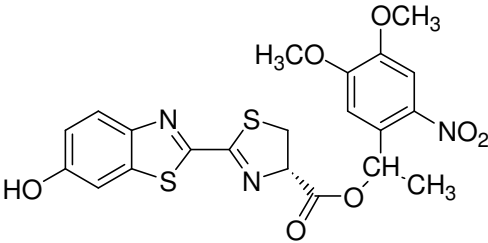


### Instruction Manual

Catalog Number	PK-CA707-10103
Description	DMNPE-caged D-luciferin is a D-luciferin ester derivative which can cross cell membranes efficiently. Once inside the cells, the ester is continuously hydrolyzed to a supply of D-luciferin. Alternatively a burst of D-luciferin is generated by UV photolysis.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} = 333 \text{ nm}$ ; Extinction coefficient: 22,000
Molecular Structure	
Molecular Weight / Molecular Formula	489.52 Da; C <sub>21</sub> H <sub>19</sub> N <sub>3</sub> O <sub>7</sub> S <sub>2</sub>
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Light yellow solid; soluble in DMF and DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	NA
References	1) <i>Antimicrobial Agents &amp; Chemotherapy</i> , 1121(Apr-2001) 2) <i>Lett Appl Microbiol</i> <b>30</b> , 223(2000) 3) <i>Biotechniques</i> <b>15</b> , 848(1993)
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.