

Product Information

Description:	Substrate for ceramide kinase (CERK), optimized for use with β -Octylglucoside mixed micellar assay
Format:	A 2-component set containing sufficient reagents for 150 \times 50 μ l assays: <ul style="list-style-type: none">• Ceramide (0.3 ml @1.25 mM)• Ceramide diluent buffer (2 \times 1.75 ml)
Stability:	At least 6 months at -20°C.
Storage:	-20°C
Substrate Preparation:	Prepare a fresh 1:4 mixture of ceramide and ceramide buffer. Sonicate the mixture 3 times for 2 min. using a bath sonicator to make a 2.5 X solution (250 μ M ceramide). A final ceramide concentration of 100 μ M is recommended. Discard any unused substrate.
Assay Conditions:	Ceramide Kinase (BPS Bioscience, #40617) reaction was performed for 25 minutes at 37°C using 100 μ M CERK substrate. Kinase-Glo [®] (Promega) was used to detect the enzyme activity.
Applications:	Useful as a substrate for CERK activity assays.

Quality Control Data

Ceramide Kinase Specific Activity

