

Company Name:
Quote #:

SAMPLE SUBMISSION FORM

Please follow the instructions below for submitting material for testing.

- 1) Label the vial with the material identification #.
- 2) Submit the material as a 10 mM solution in DMSO (100 uL) or as a pre-weighed solid (equivalent amount to make a 10 mM stock).
- 3) Include any special instruction for sample handling or storage.
- 4) Fill out the table below and send this form along with the sample.

Material ID#	Sample form (Solid/Liquid)	10 mM in DMSO? (Y/N)	For solid samples only		Screening Conc (uM)	Special Instructions
			Amount sent	Molecular Weight		

Samples submitted for analysis should be properly packaged and shipped to the following address:

FedEx/UPS/US Mail:
 Luceome Biotechnologies
 1665 E. 18th Street, # 106
 Tucson, AZ 85719
 Ph: (520) 495-0161