

## Product Information Sheet

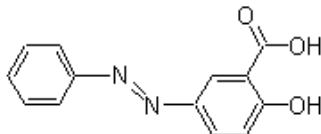
### Ordering Information

Product Number: 23054  
Product Name: Half chrysamine G  
Unit Size: 5 mg  
Storage Conditions: <-15 °C and kept from light and moisture  
Expiration Date: 12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 242.2

Chemical Structure:



Soluble in: DMSO  
Excitation Wavelength: 342 nm  
Emission Wavelength: NA

### Application Notes

A half-molecule of Chrysamine G, the well-known marker for beta-amyloid deposition, offers protection against beta-amyloid 25-35 and beta-amyloid 40-induced neuronal death at a concentration of 0.1-1  $\mu$ M. It is shown to cross the blood brain barrier. This molecule possesses minimal *in vivo* toxicity and displays low affinity to beta-amyloid40.