

## DAA

A probe for nitric oxide (NO) that forms a red ( $\lambda_{Em}$  580 nm) fluorescent precipitate after reacting with NO.



### Product attributes

CAS number	1758-68-5
Excitation/Emission	488/580 (After reaction with NO)
Molecular weight	336.32
Storage Conditions	Store at -10 to -35 °C, Protect from light, Desiccate

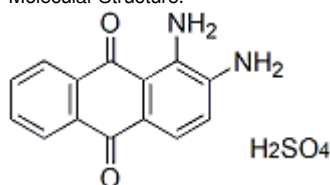
## Product Description

DAA (1,2-Diaminoanthraquinone) reacts with nitric oxide (NO) to form a red ( $\lambda_{Em}$  580 nm) fluorescent precipitate.

- Probe for nitric oxide (NO)
- Reacts with NO to form a red fluorescent precipitate
- $C_{14}H_{12}N_2O_6S$
- Yellow-orange solid soluble in DMSO

DAA has been used to detect NO in cultured neurons and in rat retinas after injury to the optic nerve.

Molecular Structure:



## References

1. Neuroreport. 9, 4051 (1998) [PMID: 9926846](#)
2. Microchim Acta 152, 35 (2005) [DOI: 10.1007/s00604-005-0420-x](#)

This datasheet was generated on January 30, 2026 at 10:20:17 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/daa-12-diaminoanthraquinone/>