



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Carboxyrhodamine 110 hydrochloride, 5 and 6 isomers



For some applications, 5-(and-6)-carboxyrhodamine (5(6)-CR 110) is a superior alternative to the widely used green fluorescent dye 5-(and-6)-carboxyfluorescein. The dye is insensitive to pH between 4 and 9 and is much more photostable than fluorescein.

Product attributes

Excitation/Emission	502/524 nm
---------------------	------------

Product Description

For some applications, 5-(and-6)-carboxyrhodamine (5(6)-CR 110) is a superior alternative to the widely used green fluorescent dye 5-(and-6)-carboxyfluorescein. The dye is insensitive to pH between 4 and 9 and is much more photostable than fluorescein.

This datasheet was generated on June 23, 2026 at 02:58:04 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-and-6-carboxyrhodamine-110-hydrochloride-56-cr-110/>