10-Acetyl-3,7-dihydroxyphenoxazine (Amplex® Red)

sicitum 680

CAS number 119171-73-2

Product attributes

Excitation/Emission 571/585 nm (end product)

Call us: 800-304-5357 Email: btinfo@biotium.com

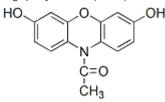
10-Acetyl-3,7-dihydroxyphenoxazine (Amplex® Red), is regarded as the best fluorogenic substrate for peroxidase and H₂O₂ because it is highly sensitive, specific and stable.

Product Description

10-Acetyl-3,7-dihydroxyphenoxazine (N-Acetyl-3,7-dihydroxyphenoxazine or ADHP), which has been marketed under the name Amplex® Red, is regarded as the best fluorogenic substrate for peroxidase because it is highly specific and stable. The substrate itself is nearly colorless and nonfluorescent until it is oxidized by H_2O_2 (reacting in a 1:1 stoichiometry) in the presence of horseradish peroxidase (HRP) to become the highly red fluorescent resorufin. H_2O_2 is produced by many enzymatic reactions, and therefore ADHP can be used to detect the activity of many different enzymes.

- Detect HRP in immunoassays
- Detect trace amounts of hydrogen peroxide (H₂O₂)
- Detect using fluorometer or microplate reader
- $\lambda_{Ex}/\lambda_{Em}$ (pH 9)= 571/585 nm (end product)
- ϵ = 54,000 (end product)
- White to off-white solid soluble in DMSO or DMF
- Store desiccated at -20°C and protect from light
- C₁₄H₁₁CNO₄
- MW: 257.24
- [119171-73-2]

High purity resorufin (10062) is also available from Biotium to be used as a standard.



Amplex Red is a registered trademark of Thermo Fisher Scientific.

References

- 1. Biosensors and Bioelectronics (2010) doi:10.1016/j.bios.2010.07.126
- 2. Cancer Biotherapy and Radiopharmaceuticals 25(4), 439 (2010).
- 3. Anal Biochem 287, 196 (2000).
- 4. J Immunol Methods 202, 133 (1997).
- 5. Anal Biochem 253, 169 (1997).

This datasheet was generated on November 17, 2025 at 11:06:41 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/10-acetyl-37-dihydroxyphenoxazine-n-acetyl-37-dihy