

DISPENSING SERVICE product release form

Requestor: NAAmp Product Manager
Catalog product

NRU: 1001106378
Internal reference: CS-ALIQ-PROD22031
Product: Red Diamond Taq®

Eurogentec Batch number:
Enzyme: 22031-P01
Reaction buffer: 22031-P01
MgCl₂: 22031-P01

Product Reference and quantity released:
TAQ-I041-1000 60 kits
TAQ-I041-5000 50 kits

Expiry date: 20/08/2026

QC results:

QC type:	Status:
Appearance (Colourless solution)	Passed / Failed
Volumique Activity ($\geq 5U/\mu l$)	Passed / Failed
Test de performance / PCR sur λ DNA	Passed / Failed
<i>q-PCR SPC</i>	<i>Supportive Data</i>
Bioburden (≤ 10 CFU/U)	Passed / Failed

- QC Comments: /
- Additional QC criteria, specific to Taq project (cf. requirements on production data - Annexe 1 of MO-DIS-005):
Passed/Failed / Not applicable*
- Approval of QC results – Pj Management:
 - Date: 26/11/2024
 - Full name + title: Axelle Lemaître – Project Leader Dx
 - Signature:



- Approval of QC results – QC Department:
 - Date: 26/11/2024
 - Full name + title: Thomas Fett – QC Assistant Manager
 - Signature:

