

DISPENSING SERVICE product release form

Requestor: NAAmp Product Manager
Catalog product

NRU: 1001106378

Internal reference: CS-ALIQ-PROD22030

Product: HGS Diamond Taq

Eurogentec Batch number:

Enzyme: 22030-P01

Reaction buffer: 22030-P01-bis

MgCl₂: 22030-P01

Product Reference and quantity released:

TAQ-I011-1000+ 138 kits

TAQ-I011-5000+ 122 kits

Expiry date: 20/03/2026

QC results:

QC type:	Status:
Appearance (Colourless solution)	Passed / Failed
Volumique Activity (≥ 5U/μl)	Passed / Failed
Test de performance / PCR sur λ DNA	Passed / Failed
q-PCR SPC	Passed / Failed
Hotstart (SYBR Green® qPCR)	Passed / Failed
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: 04/09/2024
- Full name + title: A. Lemaître – Project leader Dx
- Signature:



Approval of QC results – QC Department:

- Date: 04/09/2024
- Full name + title: T. Fett – Qc Assistant Manager
- Signature:

