

DISPENSING SERVICE product release form

Customer (Name, Institute):

Catalog product – cf Genomics Product manager

NRU: 1001072759
Product Reference and quantity delivered:

TAQ-I021-1000 150 kits
TAQ-I021-5000 100 kits
Internal reference: CS-ALIQ-PROD22028



Eurogentec Batch number: 22028-P01

Customer batch number (optional): NA

Expiry date: 08/03/2025

Standard QC results:

QC Plate Name	Negative Control IPC	Negative Control customer target	Positive Control IPC	Ct StDev
20230310- test QC Diamond - IPC	NA	NA	Passed / Failed	Passed / Failed

- QC Comments: NA
- Additional QC criteria, specific to *Taq* project (cf. requirements on production data - Annexe 1 of MO-DIS-005): Passed / Failed / Not applicable
 - Bioburden (results received on June 30th)
 - NPD report: QC (NPD)CS-ALIQ-PROD22028_Lot22028-P01.pdf on 26th of September 2023
- Approval of QC results – Pj Management:
 - Date: 09/11/2023
 - Full name + title: A. Lemaître – Project leader Oligo
 - Signature: 
- Approval of QC results – QC Department:
 - Date: 22/11/2023
 - Full name + title: N. SHARCIASSO QC MANAGER
 - Signature: 

Checklist documents:

- Protocol / product description / NA (Internal catalog product)
 - Validated by customer on: ___ / ___ / _____
 - No validation required (for recurrent orders or if it is a sample batch)
- Oligonucleotides / probes TDS – NA
- Oligonucleotides / probes QC – NA
- Mastermix Certificate of Analysis
 - Sent to customer
 - Stored on KEGT catalogue product files (for internal mastermix production)
- Other (please specify):

Shipment operator: NA