



Red Load Taq Master Lyophilisate

Lyophilized Taq Master Mix containing red gel loading dye
preloaded 8-tube strips

Cat. No.	Amount
PCR-151S-8TS	12 strips (96 reactions)
PCR-151L-8TS	60 strips (480 reactions)

For *in vitro* use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Additional Storage Conditions: Store in an aluminium-coated bag or on a dry place. Lyophilisates may hydrate at humidity levels >70 % when sealing is opened.

Shelf Life: 12 months

Form: lyophilized

Description:

Red Load Taq Master Lyophilisate is delivered in PCR reaction tube strips or 96-well plates preloaded with a complete master mix in a dry, room temperature stable format. The lyophilisate combines highest performance with convenience of use and stability. There is no need for freezing, thawing or pipetting on ice. The few remaining pipetting steps minimize the risk of errors or contaminations.

Each vial contains polymerase, dNTPs and reaction buffer with MgCl₂ required for a 20 µl PCR assay. The Red Load Taq Master Lyophilisate contains additionally an inherent red dye allowing the direct loading of the PCR reaction product onto the gel.

To perform PCR, fill up the vials with a premix of primers and PCR-grade water and add DNA template. If necessary, centrifuge to remove bubbles, vortex the vials to assure homogeneity and start cycling.,

Kit contents:

Red Load Taq Master Lyophilisate,

Preloaded lyophilisates of Taq DNA polymerase, dATP, dCTP, dGTP, dTTP, Reaction Buffer, MgCl₂, red gel loading dye and stabilizers.,

PCR grade water

Recommended 20 µl PCR assay:

forward Primer	0.2 - 1 µM
reverse Primer	0.2 - 1 µM
template DNA	1 - 50 ng
PCR-grade water	fill up to 20 µl

Recommended cycling conditions:

Initial denaturation	94 °C	2 min	1x
Denaturation	94 °C	30 sec	30x
Annealing ¹⁾	50 - 68 °C	30 sec	30x
Elongation ²⁾	72 °C	30 sec - 3 min	30x
Final elongation	72 °C	2 min	1x

¹⁾ 1) The annealing temperature depends on the melting temperature of the primers used. ,

²⁾ 2) The elongation time depends on the length of the fragments to be amplified. A time of 1 min/kb is recommended.,

**Red Load Taq Master Lyophilisate**

Lyophilized Taq Master Mix containing red gel loading dye
preloaded 8-tube strips

For optimal specificity and amplification an individual optimization of the recommended parameters may be necessary for each new template DNA and/or primer pair.