

# NZYTaq II 2× Green Master Mix

Catalogue number	Presentation
MB35801	2 x 1.25 mL (100 rxns of 50 µL)
MB35802	4 x 1.25 mL (200 rxns of 50 µL)
MB35803	20 x 1.25 mL (1000 rxns of 50 µL)
MB35805	1 x 50 mL (2000 rxns of 50 µL)

## Description

NZYTaq II 2× Green Master Mix is a premixed ready-to-use solution containing NZYTaq II DNA polymerase (NZYtech, Cat. No. MB354), which belongs to a new generation of Taq-derived DNA polymerases optimized for standard PCR applications. The master mix contains dNTPs, reaction buffer and additives at optimal concentrations and supports the robust and reliable amplification of a wide range of DNA templates up to 6 kb. MgCl<sub>2</sub> final concentration is 2.5 mM, allowing the implementation of a comprehensive range of PCR protocols. In addition, reactions assembled with NZYTaq II 2× Green Master Mix may be directly loaded onto agarose gels. The mix contains two dyes (blue and yellow) that allow monitoring the progress of the electrophoresis. NZYTaq II 2× Green Master Mix is not suitable when direct fluorescent or absorbance readings are required without prior purification of the amplified DNA from PCR. We recommend using the master mix version without dyes – NZYTaq II 2× Colourless Master Mix (NZYtech, Cat. No. MB357) – or purifying the PCR product using NZYGelpure (NZYtech, Cat. No. MB011) before performing such downstream protocols.

NZYTaq II DNA polymerase lacks 3'→5' exonuclease activity. Resulting PCR products have an A-overhang and are suitable for cloning with NZYtech's NZY-A PCR cloning kit (NZYtech, Cat. No. MB053) or NZY-A Speedy PCR cloning kit (NZYtech, Cat. No. MB137).

## Shipping & Storage Conditions

The product can be shipped from dry ice to ambient temperature. Upon arrival, all components should be stored at -85 °C to -15 °C in a constant-temperature freezer to guarantee maximal shelf life. The high thermal stability of the enzyme mixture allows it to remain stable at 4 °C or even at room-temperature for up to 4 weeks. The product will remain stable till the expiry date if stored as specified.

## Components

COMPONENT	SKU	TUBES/BOTTLES	VOLUME
NZYTaq II 2× Green Master Mix	MB35801	2	1.25 mL
	MB35802	4	1.25 mL
	MB35803	20	1.25 mL
	MB35805	1	50 mL

## Standard Protocol

### Recommendations before starting

- **Nucleic acid manipulation:** The quality and integrity of DNA templates are critical for achieving reliable PCR results. Use high-quality DNA extraction kits to isolate clean, contaminant-free DNA. To avoid degradation or contamination, handle DNA samples with care and work in designated clean areas. Consider using NZYtech Nucleases & Nucleic Acid Cleaner (Cat. No. MB48301) or DNA & RNA Cleaner (Cat. No. MB46201) to remove nucleases or residual contaminants from surfaces and materials.
- **Handling instructions:** To prevent carry-over contamination, dedicate separate areas for reaction setup, PCR amplification, and any post-PCR analysis. Never open tubes containing amplified PCR products in the PCR setup area. Always use sterile, filtered pipette tips and avoid reusing any consumables.
- **Preventing Contamination:** Implement stringent laboratory practices to avoid false positives caused by contaminants. Use sterile equipment, clean workstations regularly, and monitor assay integrity by including No-Template Controls (NTC) in each PCR run.

## Procedure

The following protocol serves as a general guideline and a starting point for any PCR procedure. Optimal reaction conditions (e.g. incubation times, temperatures and template concentration) may vary and, in particular conditions, may require further optimization. In case you need to fine-tune primer concentrations, test recommended variations provided in brackets in the table below.

1. Thaw the master mix at room temperature or on ice. Mix the master mix thoroughly by flicking the tube and inverting it. Keep the master mix on ice. Set up the PCR reaction on ice and add water first and the remaining components in the order specified in the table below. A single reaction mixture of 50  $\mu$ L should combine the following components:

	1 REACTION VOLUME
NZYTaq II 2 $\times$ Green Master Mix	25 $\mu$ L
Primers	0.25 (0.1-0.5) $\mu$ M
Template DNA	5 $\mu$ g-0.5 $\mu$ g
Nuclease-free water	up to 50 $\mu$ L
FINAL VOLUME =	50 $\mu$ L

2. Gently mix and centrifuge briefly to spin down the contents.
3. Perform PCR using the following cycling parameters:

CYCLES	TEMP.	TIME	STAGE
1	95 $^{\circ}$ C	3 min	Initial denaturation
25-35	94 $^{\circ}$ C	30 sec (¥)	Denaturation
	(*)	30 sec	Annealing
	72 $^{\circ}$ C	15-30 sec/kb (¥)	Extension
1	72 $^{\circ}$ C	5-10 min	Final Extension

(\*) Annealing temperature should be optimized for each primer set based on the primer  $T_m$ ; typically, it should be  $T_m - 5$   $^{\circ}$ C.

(¥) For DNA fragments higher than 3 kb to 6 kb in size, it may be beneficial to use 20 sec for denaturation and 30-60 sec/kb for extension.

4. Analyse the PCR products through agarose gel electrophoresis (0.7-1.2%, w/v) and visualise with GreenSafe Premium (NZYtech, Cat. No. MB132) or any other means.

## Technical Notes

**Primer design:** PCR primers generally range in length from 15–30 bases and are designed to flank the region of interest. Primers should contain 40–60% GC and care should be taken to avoid sequences that might produce internal secondary structure. The 3'-ends of the primers should not be complementary to avoid the production of primer-dimers. Primer-dimers unnecessarily delete primers from the reaction and result in an unwanted polymerase reaction that competes with the desired reaction. Avoid three G or C nucleotides in a row near the 3'-end of the primer, as this may result in non-specific primer annealing. Ideally, both primers should have nearly identical melting temperatures ( $T_m$ ), allowing their annealing with the denatured template DNA at roughly the same temperature.

**DNA template:** The optimal amount of starting material may vary depending on its quality and complexity. In general, we recommend using 10 ng to 500 ng of genomic DNA templates, although the enzyme is sensitive enough to amplify fragments from as little as 5  $\mu$ g of human gDNA, for example. Lower amounts of template may be used for amplification of less complex DNA (typically 1-20 ng). When using a cDNA synthesis reaction as template do not exceed 10% of the final PCR reaction volume.

**PCR controls:** The reliability of the data may be affected by the presence of contaminating DNA. It is strongly recommended to include a no-template control reaction in the PCR design, replacing template DNA with nuclease-free PCR-grade water. Additionally, include a positive control to serve as a reference for ensuring the correct functioning of the PCR reaction. The positive control should exhibit the expected amplification profile, confirming the assay's ability to accurately amplify the target sequence.

**Intercalating Dyes:** The master mix contains both a blue and a yellow dye. The blue dye migrates at the same rate as a 3–5 kb DNA fragment in a 1% (w/v) agarose gel. The yellow dye migrates faster than primers (<50 bp) in a 1% (w/v) agarose gel. This mix is not suitable for direct fluorescence or absorbance measurements without prior purification of the amplified DNA from the PCR.

## Quality control assays

### Genomic DNA contamination

The product must be free of any detectable DNA contamination as evaluated through PCR.

## Nuclease assays

To test for DNase contamination, 0.2-0.3 µg of pNZY28 plasmid DNA are incubated with the master mix for 14-16 h at 37 °C. Following incubation, the nucleic acids are visualized on a GreenSafe-stained agarose gel. There must be no visible nicking or cutting of the nucleic acids.

## Functional assay

NZYTaq II 2× Green Master Mix is tested for performance in a polymerase chain reaction (PCR) for the amplification of different-sized DNA fragments (1 and 5 kb) from human genomic DNA. The resulting PCR products are visualized as single bands in a GreenSafe Premium-stained agarose gel.

## Troubleshooting

Troubleshooting is often a systematic, meticulous process where varying one parameter at a time and evaluating impacts can unveil the root cause of issues. These adjusted suggestions, incorporating a blend of specificity and exploratory approaches, aim to enhance the clarity and actionability of your troubleshooting guide. Should any other technical or procedural aspects require attention, your feedback and additional information will always be welcomed.

<b>NO PRODUCT AMPLIFICATION OR LOW YIELD</b>
<ul style="list-style-type: none"><li>• <b>Inadequate annealing temperature</b></li></ul>
The reaction mix composition may affect the melting properties of primers and DNA. Adjust the annealing temperature to accommodate the primer with the lowest melting temperature (5 ° to 10 °C lower than T <sub>m</sub> ).
<ul style="list-style-type: none"><li>• <b>Presence of PCR inhibitors</b></li></ul>
Some DNA isolation procedures, particularly genomic DNA isolation, can result in the co-purification of PCR inhibitors. Reduce the volume of template DNA in reaction or dilute template DNA prior to adding to the reaction. Diluting samples even 1:10,000 has been shown to be effective in improving results, depending on initial DNA concentration.
<ul style="list-style-type: none"><li>• <b>Template DNA damaged or degraded</b></li></ul>
An intact, high-quality template is essential to achieve a reliable amplification, especially from large DNA fragments. Extreme care must be taken in the preparation and handling of DNA. Always use purified high-quality DNA as template.
<ul style="list-style-type: none"><li>• <b>Contamination with DNases</b></li></ul>
Ensure that all labware, including pipettes, tubes, and containers, is clean and free from residual DNase contamination. Disinfect laboratory surfaces with the RNase & DNase Cleaner (NZYtech, Cat. No. MB463). Use autoclaved or sterile equipment whenever possible. Use DNase-free water. Change gloves frequently.
<ul style="list-style-type: none"><li>• <b>Concentration of Mg<sup>2+</sup> is too low</b></li></ul>
Mg <sup>2+</sup> is included in the Master Mix at a final concentration of 2.5 mM, which is sufficient for most targets. For some targets, higher Mg <sup>2+</sup> concentration may be required. Titrate from 2.5 mM to 4 mM (final concentration) in 0.5 mM increments. (Note: MgCl <sub>2</sub> is not provided in separate tubes).
<b>PRESENCE OF NON-SPECIFIC BANDS</b>
<ul style="list-style-type: none"><li>• <b>Non-specific annealing of primers</b></li></ul>
Adjust annealing conditions and/or design another set of primers, by increasing the length and avoiding complementary sequences.
<ul style="list-style-type: none"><li>• <b>Primer degradation</b></li></ul>
Check the quality and concentration of primer solutions. We recommend preparing small volume working aliquots from the stock solution. Avoid using primers subjected to multiple freezing-thawing cycles.

*This master mix is manufactured under stringent quality standards and complies with ISO 9001 and ISO 13485 certifications for research and diagnostic-grade reagents.*

For life science research only. Not for use in diagnostic procedures.