

Mixing of Highly Viscous Liquids Using Eppendorf MixMate®

Wolf Wente¹, Daniel Dauch² and Lars Borrmann²

¹Eppendorf Instrumente GmbH, Hamburg, ²Eppendorf SE, Hamburg, Germany

Abstract

Viscous liquids such as glycerol or polyethylene glycol (PEG) are often used as a nonaqueous solvent for drugs, vaccines or dyes or serve as antifreeze and stabilizing agents in protein solutions. Because of their high viscosity these solvents are hard to mix and place specific demands on the performance of laboratory mixers.

The aim of this work was to analyze the performance of Eppendorf benchtop mixer MixMate in mixing highly viscous solutions. Therefore, we determined experimentally the maximum glycerol concentration being efficiently mixed by MixMate in different plate and tube formats. Results presented herein show that MixMate is able to efficiently mix glycerol solutions up to 90 % at 20 °C which corresponds to a viscosity of 219 Centipoise (cP), i.e., approx. 219-fold the viscosity of water. The best mixing performance was found for vessels with a



wide diameter and flat bottom. In general, the mixing performance of MixMate allows fast and efficient mixing of aqueous solutions as well as highly viscous samples.

Introduction

MixMate is a high speed mixer that was especially developed for fast and controlled mixing of small reaction volumes in all types of plates, including 384-well plates.

As proven in previous publications, MixMate mixes all kinds of molecular biological samples within 1 minute, independent of sample properties or plate geometry [1]. While most samples and reaction buffers used in molecular biology are based on aqueous solutions having a viscosity

comparable to water we also wanted to test the capability of MixMate in mixing highly viscous solutions. These highly viscous solutions such as glycerol, propylene glycol or polyethylene glycol (PEG) are often used in pharmaceutical or food industry as nonaqueous solvents for drugs, vaccines or dyes. In addition, with molecular biology enzyme solutions, these substances serve as antifreeze and stabilizing agents.

Viscosity, commonly perceived as “thickness”, is a measure that describes the resistance of a fluid to deform under shear stress. The viscosity varies with the substance, its concentration and with temperature. The higher the temperature, the lower the viscosity.

There are different physical units in use to express the dynamic viscosity of a solution: The International System of Units (SI unit) expresses viscosity in pascal-second (Pa·s) whereas the centimeter-gram-second system (CGS) uses poise (P), or more commonly centipoise (cP). Centipoise is commonly used because water has a viscosity of 1.0020 cP (at 20 °C) [2,3]. The relation between both units is:

$$1 \text{ Pa}\cdot\text{s} = 10 \text{ P}$$

$$1 \text{ mPa}\cdot\text{s} = 1 \text{ cP}$$

Materials and Methods

The mixing performance of MixMate was tested with different vessel types (micro test tubes and 96-well DWP, MTP and PCR plates; please refer to Table 1 for details) and different filling volumes (12.5, 25, 50 and 75 % of the recommended working volume).

For the assay, glycerol (Rotipuran® ≥ 99.5 %, p.a. water-free, Carl Roth GmbH, Karlsruhe, Germany) solutions of different concentrations (50, 60, 70, 75, 80, 85, 90, 95 % (v/v) in dH₂O) were pipetted to the bottom of the vessels and were carefully overlaid with 1/30 volume of Ponceau-4R colored dH₂O (Figure 1). After 1 min mixing with Eppendorf MixMate at maximum speed (3,000 rpm for MTP and PCR plates; 2,000 rpm for tubes and DWP) the vessels were visually checked for completely mixed glycerol solutions. A solution was rated as completely mixed if the solution was stained homogeneously red without any cords (Figure 1).

All experiments were performed at 20 °C room temperature. Solutions, plates, tubes, MixMate, pipettes and tips were equilibrated to room temperature.

For example, the viscosity of some undiluted substances are:

> Glycerol: 1410 mPa·s at 20 °C

> PEG-200 (MW 190 – 210): 50 mPa·s at 25 °C

> PEG-300 (MW 280 – 320): 70 mPa·s at 25 °C

> PEG-400 (MW 380 – 420): 90 mPa·s at 25 °C

> PEG-600 (MW 570 – 630): 135 mPa·s at 25 °C

> Propylene glycol: 40.4 mPa·s at 25 °C

> Triton® X-100: 240 mPa·s at 25 °C

In this work we analyzed the performance of MixMate in mixing highly viscous solutions. Therefore, we determined experimentally the maximum glycerol concentration being efficiently mixed by Eppendorf MixMate in different kind of tubes and 96-well plates and at different filling volumes. All experiments were performed at 20 °C. The tested glycerol concentrations were afterwards correlated to their corresponding viscosity in mPa·s as given within the literature [4].

Results and Discussion

Previous publications have shown that the unique 2DMix-Control technology of MixMate enables fast mixing of small reaction volumes independent of sample properties or vessel geometry [1]. Even difficult mixing applications such as resuspension of bacterial cell pellets can be achieved with MixMate in less than 1 minute [6]. Furthermore, the 2Dmixing stroke ensures a controlled mixing motion which reduces the risk for lid-wetting or cross-contamination [7].



Figure 1: Experimental setup before (left) and after (right) mixing with MixMate. Left: Glycerol solutions overlaid with 1/30 volume of red colored water. Right: A completely mixed solution as indicated by the homogeneously red staining without any cords.

In this Application Note we analyzed the mixing performance of MixMate by mixing glycerol solutions in different kind of tubes and 96-well plates. The experiments showed that MixMate can efficiently mix glycerol solution up to 90 % in less than 1 minute (Table 1). According to data provided in the literature a 90 % glycerol concentration corresponds to a viscosity of 219 mPa.s, 219-fold the viscosity of water [4,5].

As can be seen by the differences between mixing results in 1.5 ml and 2.0 ml tubes and micro test plates with round and flat bottom, the shape of the vessel bottom largely influences the mixing performance. Flat bottom vessels

were easier to mix than those with round bottom or conical shape.

Looking at the relationship between sample viscosity and total sample volume, the data show that the best performance at high viscosities was achieved with Eppendorf Plates Deepwell 96 2 ml (DWP 96/2000) as up to 500 μ l of 90 % glycerol solution (viscosity = 219 mPa.s) were efficiently mixed by MixMate. These data prove that MixMate, besides its superior performance in mixing aqueous solutions, also enables efficient mixing of highly viscous samples.

| Tubes | | Plates (96-well) | | | | |
|-----------------------|--------------------------------------|------------------------------------------|--------------------------------------------|---------------------------------------|--------------------------------------|-------------------------------------------|
| | Micro test tubes (max. 2,000 rpm) | Deepwell plate (DWP) (max. 2,000 rpm) | Micro test plate (MTP) (max. 3,000 rpm) | PCR plate (max. 3,000 rpm) | | |
| | | | | | | |
| | | | | | | |
| Total volume: | Eppendorf Tube Safe-Lock 2.0 ml | Eppendorf Plate Deepwell 96/2000 µl | Eppendorf Plate Deepwell 96/1000 µl | Nunc F96 MicroWell™ Plates (# 260836) | Nunc 96 MicroWell™ Plates (# 249570) | Eppendorf twin.tec® PCR Plate 96, skirted |
| Working volume: | 2,000 µl | 2,400 µl | 1,200 µl | 400 µl | 300 µl | 150 µl |
| Bottom shape: | 2,000 µl | 2,000 µl | 1,000 µl | 200 µl | 150 µl | 150 µl |
| 12.5 % filling volume | round | RecoverMax conical | RecoverMax conical | flat | U-bottom | conical |
| 25 % filling volume | 85 % (109 mPa s) | 90 % (219 mPa s) | 85 % (109 mPa s) | 90 % (219 mPa s) | 90 % (219 mPa s) | 80 % (60 mPa s) |
| 50 % filling volume | 85 % (109 mPa s) | 90 % (219 mPa s) | 80 % (60 mPa s) | 90 % (219 mPa s) | 90 % (219 mPa s) | 80 % (60 mPa s) |
| 75 % filling volume | 80 % (60 mPa s) | 80 % (60 mPa s) | 70 % (22 mPa s) | 90 % (219 mPa s) | 85 % (109 mPa s) | 80 % (60 mPa s) |
| | 75 % (35 mPa s) | 70 % (22 mPa s) | 50 % (6 mPa s) | 90 % (219 mPa s) | 85 % (109 mPa s) | 50 % (6 mPa s) |

Table 1: Mixing of highly viscous solutions in tubes and 96-well plates using Eppendorf MixMate. Vessels were filled with glycerol solution up to 12.5, 25, 50 and 75 % of the recommended working volume. The maximum glycerol concentration (in percent) that was completely mixed after 1 minute at either 2,000 rpm (tubes and DWP) or 3,000 rpm (MTP and PCR plate) is listed. In parentheses the viscosity of the aqueous glycerol solutions in millipascal-seconds (mPa s) is given. All experiments were performed at a temperature of 20°C.

References

- [1] Osterhoff C, Dauch D, Mueller P, Borrmann L. Comparison of mixing performance in 96- and 384-well plates of Eppendorf MixMate and competitor devices. Eppendorf Application Note 130, 2006.
- [2] Swindells JF, Coe JR, Godfrey TB. Absolute viscosity of water at 20 °C. J. Res. Nat. Bur. Stand. 1952; 48(1):1-31.
- [3] Roscoe R and Bainbridge W. Viscosity Determination by the Oscillating Vessel Method II: The Viscosity of Water at 20 °C. Proc. Phys. Soc. 1958; 72:585-595.
- [4] Segur JB and Oberstar HE. Viscosity of Glycerol and Its Aqueous Solutions. Industrial and Engineering Chemistry. 1951; 43(9):2117 – 2120.
- [5] The Dow Chemical Company (www.dow.com/glycerine). Brochure: Discover the origins of some of the world's most consistently pure products: Synthetic glycerine products. Form No. 115-00645-799 SMG, July 1999
- [6] Osterhoff C, Mueller P, Borrmann L. Eppendorf MixMate – Resuspension of bacteria pellets in deepwell plates (96- and 384-well) and micro test tubes. Eppendorf Application Note 131, 2006.
- [7] Osterhoff C, Mueller P, Borrmann L. Eppendorf MixMate – Experimental evidence of controlled mixing using a PCR-based chessboard assay. Eppendorf Application Note 129, 2006.

Ordering information

| Description | Order no. International | Order no. North America |
|--------------------------------------------------------------------------------|-----------------------------|-------------------------|
| MixMate® , basic device without tube holders | | |
| 230 V, 50–60 Hz | 5353 000 510 | 2231000804 |
| 110 V, 50–60 Hz | 5353 000 529 | |
| Eppendorf Safe-Lock Tubes 1.5 ml , Eppendorf quality, colorless | 0030 120 086 (1,000 pcs) | 022363204 (500 pcs) |
| Eppendorf Safe-Lock Tubes 2.0 ml , Eppendorf quality, colorless | 0030 120 094 (1,000 pcs) | 022363352 (500 pcs) |
| Deepwell Plate 96/1000 µL , PCR clean, Blue, 20 plates (5 bags of 4) | 0030 501 241 | 951032689 |
| Deepwell Plate 96/2000 µL , PCR clean, Blue, 20 plates (5 bags of 4) | 0030 501 349 | 951033481 |
| Eppendorf twin.tec® PCR Plate 96, skirted 150 µL , colorless, 25 pcs | 0030 128 508 | 951020401 |

Your local distributor: www.eppendorf.com/contact
 Eppendorf SE · 22339 · Hamburg · Germany
eppendorf@eppendorf.com · www.eppendorf.com

www.eppendorf.com

Rotipuran® is a registered trademark of Carl Roth GmbH, Karlsruhe, Germany.

Triton® is a registered trademark of Union Carbide Co.

Eppendorf®, the Eppendorf Brand Design, MixMate®, Mastercycler® and Eppendorf twin.tec® are registered trademarks of Eppendorf SE, Germany. All rights reserved, including graphics and images. Eppendorf SE reserves the right to modify its products and services at any time. This white paper is subject to change without notice. Although prepared to ensure accuracy, Eppendorf SE assumes no liability for errors, or for any damages resulting from the application or use of this information. Reading this white paper alone cannot as such provide for or replace reading and respecting the current version of the operating manual.

All rights reserved, including graphics and images. Copyright © 2024 by Eppendorf SE, Germany. Order No: AA01 038 720/EN2