

HAAKE CaBER 1 – binary polymer mixtures

Polystyrene, blends of standards,

$M_w = 1.8 \text{ Mio g/mol} + M_w = 13 \text{ Mio g/mol}$

Rheology Application Notes

CaBER test result

At a constant concentration of 250 ppm in Boger fluids binary mixtures can be differentiated in extensional flow.

In shear flow differences between in mixing ratio cannot be detected.

Conclusion

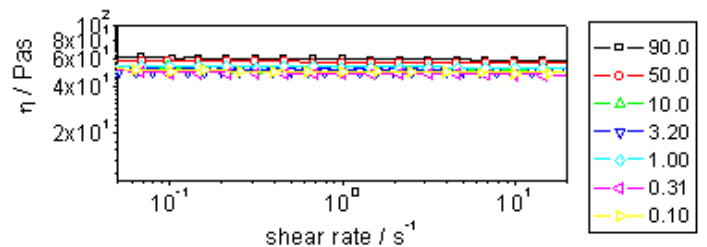
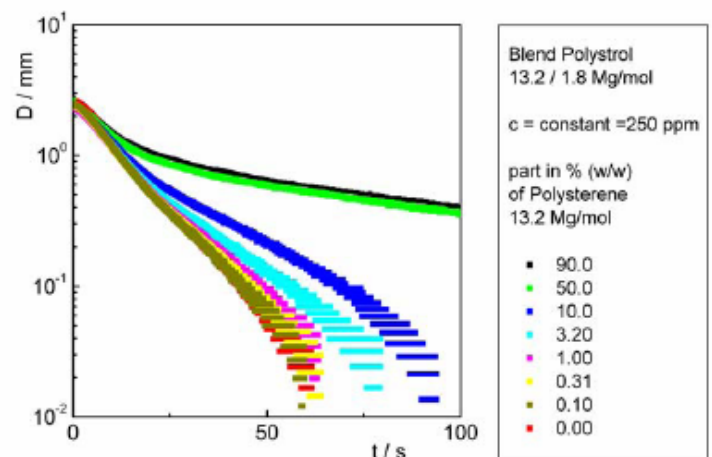
The influence of molecular weight of polymers is more sensitive in the CaBER measurement.

Practical applications

Industrial processes where small changes in polymerization process or recipe may influence application behaviour.

References

Clasen/ University of Hamburg
RheoFuture 2004



**Thermo Fisher Scientific
Process Instruments**

International/Germany
Dieselstr. 4,
76227 Karlsruhe
Tel. +49(0)721 40 94-444
info.mc.de@thermofisher.com

Benelux
Tel. +31 (0) 76 5 87 98 88
info.mc.nl@thermofisher.com

China
Tel. +86 (21) 68 65 45 88
info.china@thermofisher.com

France
Tel. +33 (0) 1 60 92 48 00
info.mc.fr@thermofisher.com

India
Tel. +91 (22) 27 78 11 01
info.mc.in@thermofisher.com

United Kingdom
Tel. +44 (0) 1785 81 36 48
info.mc.uk@thermofisher.com

USA
Tel. 603 436 9444
info.mc.us@thermofisher.com

www.thermo.com/mc

V214_07.04.

© 2004/07 Thermo Fisher Scientific -
All rights reserved - This document is
for informational purposes only and is
subject to change without notice.