

Comment on the AGM from the Board of Directors regarding item 4.1.c (re-election of Heinrich Fischer as member of the Board of Directors).

The Board of Directors informs that, while Heinrich Fischer is proposed to be re-elected as a member of the Board, he is not the designated Chairman of the Audit Committee for the upcoming term 2019/2020. The Board will make a change to the Chair of the Audit Committee after the AGM in April 2019 to one of the other Board Members.

The Chairman of the Board

Maennedorf, 28th March 2019

