

Official Abcam poster competition terms and conditions

Effective November 2017

1. The contest is open to persons aged 21 years or over, except employees of, and consultants to Abcam, their families, government employees and any third party associated with administration of the contest.
2. The contest is free to enter and no purchase is necessary. VOID WHERE PROHIBITED BY LAW.
3. All entries must be submitted and presented at the Adult Neurogenesis 2018 held at the Center for Regenerative Therapies Dresden/TU Dresden on May 2-4, 2018. Entries received after this time will not be accepted.
4. The winner will be selected by a judging panel and announced at the end of the meeting. The winners will receive Amazon gift vouchers as the prize, to the amounts of 100 EUR, 100 GBP or 100 USD – first prize; 50 EUR, 50 GBP, or 50 USD – second prize; 25 EUR, 25 GBP or 25 USD – third prize.
5. The prize is non-exchangeable, non-transferable and no cash alternative is offered.
6. The decision of the judges regarding any aspect of the contest is final and binding and no correspondence will be entered into about it.
7. By entering this competition, you are agreeing to be bound by these terms and conditions, and you are representing and warranting to Abcam that you meet the eligibility requirements of the contest. You are also agreeing to be contacted by Abcam in relation to the competition. We will not share your details with anyone else except the winner's announcement.
8. By entering this competition, you agree to receive further information regarding Abcam products and services. You can opt out of receiving communications from us at any time by unsubscribing from our newsletter.
9. Insofar as is permitted by law, Abcam reserves the right to hold void, cancel, suspend, or amend the contest where it becomes necessary to do so.
10. Insofar as is permitted by law, Abcam, its agents or distributors will not in any circumstances be responsible or liable to compensate the winner or accept any liability for any loss, damage, personal injury or death occurring as a result of taking up the prize except where it is caused by the negligence of Abcam, its agents or distributors or that of their employees. Your statutory rights are not affected.
11. Personal data supplied during this contest maybe passed onto third party suppliers only insofar as required for fulfilment/delivery/arrangement of the prize.
12. The voucher credit must be used within the expiry date listed on the voucher
13. This offer is void if it conflicts with the policy of the recipient.
14. The contest will be governed by law of England and Wales and entrants to the contest agree to submit to binding arbitration to resolve any disputes relating to the contest.
15. Abcam plc located at 330 Cambridge Science Park, Cambridge CB4 0FL is the sponsor of this contest.