

## TERMS AND CONDITIONS OF MALVERN PANALYTICAL SCIENTIFIC AWARD 2022

### **1. Prize designation**

The prize is awarded for outstanding research, published by scientists at an early stage of their research career.

### **2. Eligibility of entrants**

Entrants must be the main contributor (the first author) to the publication and research in order to be eligible for receiving the Malvern Panalytical Scientific Award. An entrant should be less than 35 years old at the moment of application submission.

Publications with authors or co-authors that are Malvern Panalytical employees or any of its subsidiaries employees or their family or household members are not eligible for the award. No persons in the employ of a company that manufactures any product or solution that is in competition with those marketed by Malvern Panalytical are eligible to apply.

### **3. Basis for the award**

The award is made for research excellence in a field where a significant thrust of the investigation that is described in a publication that required the use of one or more products or solutions produced or marketed by Malvern Panalytical (<https://www.malvernpanalytical.com/en/products#tab-3>)

Suitable research areas include, for example:

- i) Structural characterization and structure-property relationship
- ii) Characterization of materials and processes that are relevant to energy storage and energy conversion (i.e. novel battery materials, fuel cells, super capacitors, etc.)
- iii) Characterization of thin films and multilayer systems
- iv) Characterization of semiconductor and electronics related materials and processes
- v) Characterization of metal alloys and powders and/or additive manufacturing processes
- vi) Characterization of nanomaterials and processes for preparing them
- vii) Characterization of organic and inorganic chemicals (i.e. polymers, MOFs, composites, catalysts, pharmaceuticals....)
- viii) Biologics: protein-based therapeutics, vaccines, cell and gene therapies, viral and non-viral vectors
- ix) Materials characterization, experimental and industrial methods in geosciences and mining
- x) Characterization materials and methods related to environmental protection

### **4. Recipient of the award**

The award is made to the primary (the first) author of a publication that:

- i) is published in a journal recognized by the Institute for Scientific Information (ISI) citation index and where the journal has an impact factor (IF) greater than 2.5 That is to say in an ISI recognized journal where the average number of citations per article is greater than 2.5 at the time that applications close on 30th September 2022.

ii) has been published in print (i.e. hardcopy) between 1st July 2020 and 31st December 2021, or officially accepted for publication.

By 'primary author' we mean the first of the names listed as authors on a publication. If the entrant is not the primary author or is not listed, the entrant must provide an explanation on why the entrant is not listed as the primary author, e.g. that the authors are listed in alphabetic order or that the first two authors have contributed equally to the work.

Applications that have been recognized by a similar award elsewhere are still eligible for the Malvern Panalytical Award.

### **5. Application for the award**

Entrants must submit their own application. Facilities for the submission of applications will be made available by Malvern Panalytical B.V.

Entrants must submit one paper covering the details of their methods and findings.

Articles on different research projects by the same primary author should be submitted on separate forms and will be counted as separate applications.

The submissions should be written in the English language. In cases where the original publication was not in English, an English translation should accompany the submission. The translation should include figure captions, legends, footnotes and supporting information. The award will be based on the English version and as such the onus for correct translation rests on the entrant.

Along with the submission of the journal article for evaluation, each application should also specify:

- i) Current contact details of the primary author
- ii) A comprehensive description of the experimental setup and instrumentation that was used
- iii) An undertaking that every other author listed on the article agrees that:
  - a. the article has been submitted for the award with their consent
  - b. the article exists in the public domain or proper permissions have been obtained from the article publisher if required so.
  - c. Malvern Panalytical will not be held responsible for any violation of copyright, patent, intellectual property, trade secret or any third-party rights
  - d. Malvern Panalytical may publicize the submissions that were short listed in the selection process and the award winner and any details made with the submission
  - e. contact details that are provided may be used as avenues to provide information of Malvern Panalytical products and solutions

### **6. Selection process**

All applications that are submitted will undergo an initial sorting and vetting process by an internal Malvern Panalytical committee and their number will be reduced to the 10 most worthy applications. The initial vetting process will seek to establish the eligibility of the application based on the criteria set out in this document. To further reduce the entries, the committee may use whatever criteria they deem fit, while placing the priority on the highest scientific value.

The committee will consist of scientists from within Malvern Panalytical or appointed by Malvern Panalytical.

The selection committee will in its sole discretion select the best entry out of above mentioned 10 entries. Preference will be given to entries:

- i) Of high scientific merit

ii) That show innovative thinking or a novel approach

Malvern Panalytical's decision is final in all matters relating to the prize draw and no correspondence or discussion will be entered into regarding the outcome. Malvern Panalytical reserves the right to check the validity of any entry and the identity of any entrant or winner and to withhold the award if Malvern Panalytical reasonably considers the entry to be in breach of these prize draw rules.

#### **8. Components and amount of the prize**

The prize consists of a cheque for the sum of € 5,000. In each case the prize money is paid to the primary author of the article, and the first author confirms that there will be no any claims made by other authors or any other party.

The tax responsibilities lie with the recipient of the award and the appropriate tax is to be paid from the € 5,000 award with no tax provision paid over and above this amount by Malvern Panalytical.

#### **9. Privacy policy**

Any data relating to an entrant received as part of the prize will be processed in accordance with Malvern Panalytical's privacy policy and data protection legislation currently in force in the Netherlands.

#### **10. Winning the prize**

Entrants accept that Malvern Panalytical will have the right, without additional permission, to use the winners' names and details of the country in which they reside for the purposes of announcing the 10 short-listed submissions and the winner of this prize. Entrants also agree to make themselves available should they be short listed or the winner, to receive the prize and present their research to an audience constituted by Malvern Panalytical.

#### **11. Registered office**

The promoter of the Malvern Panalytical Award is Malvern Panalytical B.V. whose registered office is at Lelyweg 1, Almelo7602EA, Netherlands.

#### **12. Entries**

Malvern Panalytical cannot be held responsible for a breakdown in communications systems or networks that may result in an entry not being registered, accepted or processed or if entries are incomplete, delayed, damaged, wrongly delivered or not received for any reason.

#### **13. Modification of rules**

In the event of unforeseen circumstances, Malvern Panalytical reserves the right to modify these rules prior to the closing date for entries on September 30th, 2022.

#### **14. Acceptance of rules**

By entering the prize draw, entrants will be deemed to be bound by and have accepted these rules.

#### **15. Legal basis**

The prize draw and these rules will be governed by the laws of the Netherlands and any disputes will be subject to the non-exclusive jurisdiction of the courts of the Netherlands.