

ANNEX 1 to the Sponsoring Agreement Beckman - iGEM

OFFICIAL RULES FOR THE APPLICATION FOR A BECKMAN COULTER GRANT FOR THE PARTICIPATING TEAMS:

Eligibility

The applicant must submit the iGEM Grant application (the "Grant") via the online link no later than June 12th, 2021. All applications must be submitted in the English language. The Grant is limited to approved iGEM teams, applicants who are at least 18 years of age or older, and who have submitted a complete grant application. This Grant program is void where prohibited by law. Individuals or entities classified as healthcare professionals or healthcare organizations under any applicable laws or industry codes of ethics may not participate in this program as this program is exclusively aiming at research use only projects. In the application the applicant must name the institution (the "Institution") in which lab/premises the applicant's team will conduct and execute the project. The applicant will also name the mentor, who will, on behalf of the Institution, be mentoring the project for the applicant's team.

Beckman Coulter will provide the Grant only to the winning applicant team. The person submitting the application (the "Applicant") is entering on behalf of their team, which will be the recipient of the Grant (the "Grant Recipient"). By submitting an application for a Grant, the Applicant certifies that he or she has obtained any and all permissions required from the Institution to conduct and execute the project in their lab/premises, including to have the grant being shipped to the Institution, and is not violating any policies of the Institution in doing so.

Prizes

All information about how to submit or participate to obtain any of the Beckman Coulter prizes is indicated in our iGEM Beckman Coulter webpage.

In essence, all iGEM teams can subscribe to a general Beckman products with Q&A presentation by online registration and will have free access to the recording.

Amongst all applicants the following prizes will be drawn:

- **Two iGEM teams will be awarded a 500 USD grant voucher if they answer correctly Microbioreactor questions and after lottery draw of the correct answers.** Teams can use the vouchers to order products from our genomics catalog. Upon ordering Teams will need to justify the use of these products in their project. Valuation is based on list price. Reagents will be sent to the Institution named by the applicant.
- **Six iGEM teams will be awarded a Genomics reagent Grant in the value of 2,500 USD (or local currency equivalent) in free genomics catalog products and 3 hours of application support from Beckman Coulter.** Valuation is based on list price. Reagents will be sent to the Institution named by the applicant.
- **iGEM teams (up to 6) will be awarded a Genomics device "EMnetik" as a loan** for the duration of the competition. It must be returned after the competition no later than January 2022.
- **One iGEM team will have access to the use of an Echo liquid Handler, either in presence or as a distance virtual activity by the help of one trained technician. This technical assistance will be limited to 5 days of operational use of the Echo.**

ANNEX 1 to the Sponsoring Agreement Beckman - iGEM

•

Grant selection will be done according to the project relevant or proven need to use the indicated genomics reagent or device. Applications will be received until June 12th 2021 at 11:59 PM EDT and awardees will be notified on Monday June 28th by email.

The teams must choose the reagents from our genomics catalog. The grant is valid for multiple uses with the total value not to exceed \$2,500 (USD) or currency equivalent. Shipping and handling will be managed by Beckman Coulter Field sales department and coordinated by Field marketing. The EMnetik device shipping will be handled by our marketing development scientists.

The Grant award must be redeemed by November 31, 2021. All taxes, including, without limitation, federal, state, and local taxes on or in connection with the Grant award are the sole responsibility of the Grant Recipient. No cash award alternative to the Grant will be awarded and the Grant is not transferable. Beckman reserves the right to substitute any award of equal value in its sole discretion.

Selection of Grant Recipients

One entry per Grant Recipient.

The Grant awards will be chosen by Beckman Coulter scientific personnel, in their sole discretion, and will depend on the following criteria:

- Best fit of the project to the use of the requested material
- Scientific relevance and level of innovation of the project
- Feasibility of the project within the competition stated period

In the event the Grant Recipient cannot receive the Grant award for any reason, the Grant award shall be awarded to an alternate Grant Recipient.

However, Beckman Coulter will have no influence in the decision of the iGEM on what team will be winning the competition.

Data Processing

To execute the awards, Beckman Coulter will collect certain personal data of the applicants. The purpose of processing such personal data is solely selecting, rewarding and informing the winners of the Grant award. Personal data processed includes the following data of the Applicants: first name, last name, and email address. The data is collected from the information obtained from the Applicant. There are no other sources of information. The data file can only be accessed by Beckman Coulter and their employees and representatives to the extent of their duties. Personal data may be disclosed and transferred outside the European Union (EU)/European Economic Area (EEA) for the purpose of processing personal data disclosed by the Applicant. If personal data is transferred outside the EU/EEA, the Sponsor will ensure that appropriate contractual measures (e.g. using standard contractual clauses of the EU Commission) are applied to the transfer, the transfer is lawful and that the processing and confidentiality are in compliance with the applicable laws. The data is retained for as long as it is necessary for choosing the Grant Recipients, informing the Grant Recipients and delivering of the Grant award. The data will be discarded directly afterward.

Applicants consent to their identity being revealed in announcements via Beckman Coulter's internal communications channels. For further information, please refer to Beckman Coulter's privacy policy available at <https://www.beckman.com/about-us/global-privacy-notice>.

ANNEX 1 to the Sponsoring Agreement Beckman - iGEM

Other Terms and Conditions

Beckman Coulter reserves the right to cancel, suspend and/or modify the Grant program in its sole discretion. By participating in the Grant program, Applicants agree to release and hold harmless Beckman Coulter and its affiliates and subsidiaries, and their employees, officers, directors, agents, and representatives from and against any claim or cause of action arising out of participation in the Grant program, including any errors in the administration of the Grant, printing errors, and technical errors, as well from receipt or use of any Grant awarded and the disclosure of their respective names in connection with Beckman Coulter's promotion of the Grant program.

Return of any Grant award notification as undeliverable will result in forfeiture of the Grant award. Legal recourse is excluded. Beckman Coulter and its affiliates and subsidiaries, and their officers, directors, employees, agents and representatives are not responsible and shall not be liable for: (a) electronic, hardware or software program, network, internet or computer malfunctions, failures or difficulties of any kind, including without limitation server malfunctions or by any error (human or otherwise) which may occur in the processing of Grant applications; or (b) any condition caused by events beyond the control of the Sponsor that may cause the application and award process to be disrupted or corrupted. Any attempt by an Applicant to deliberately damage any website or undermine the legitimate operation of the Grant program may be a violation of criminal and civil laws. Should such an attempt be made, the Sponsor reserves the right to seek damages from such person to the extent permitted by law.