



BECKMAN COULTER CARES AWARD



WE CARE THAT YOU CARE

KEY DATES



NOMINATIONS

1 DEC | 2016 -
14 APRIL | 2017

Call for nominations
window



COMMUNITY VOTING

14 APRIL - 31 AUG | 2017

Community voting
window



SHORT LISTING

31 AUG - 9 NOV | 2017

International Panel
of Judges reviews
candidates and
shortlists



AWARD RECIPIENT

1 DEC | 2017

Beckman Coulter
announce the 2017
Award recipient

As the nomination process requires a significant commitment of time and energy on your part, we encourage you to review the guidelines and forms carefully before proceeding

The annual cycle of The Beckman Coulter CARES Award begins in the World AIDS Day 2016, 1st December with a 'Call for Nominations'. Nominees are brought to the attention of the Award by a nominating team consisting of a Nominee and a Nominator. Nominees cannot self-nominate. The Nominator acts as the point of contact between the Award and the nominee during the nomination and evaluation process.

The 'Call' opens December 1st 2016 and closes 14th April 2017 to allow nominating teams to adequately prepare their nominations. As the nomination process requires a significant commitment of time and energy on your part, we encourage teams to review the **nomination rules** and the **nomination forms** carefully before proceeding. Please refer to the **FAQs** page with questions and concerns and how to deliver your nomination for delivery instructions.

Between April 15th and August 31st 2017 Nominees are then listed on the CARES Award website to enable voting by the HIV /AIDS community.

The CARES Award International Panel of Judges then shortlists and chooses a winner between August 31st - November 9th 2017.

The Award winner will be announced at the International Conference on AIDS and STIs in Africa (ICASA) which will be held in Abidjan, Cote D'Ivoire, 4th - 9th December, 2017

All nominations will be duly acknowledged by email, to all members of the nominating team, with a copy to the Nominee.