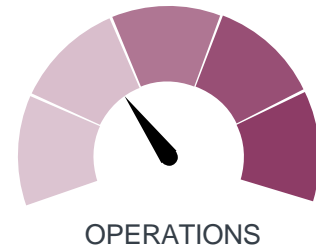
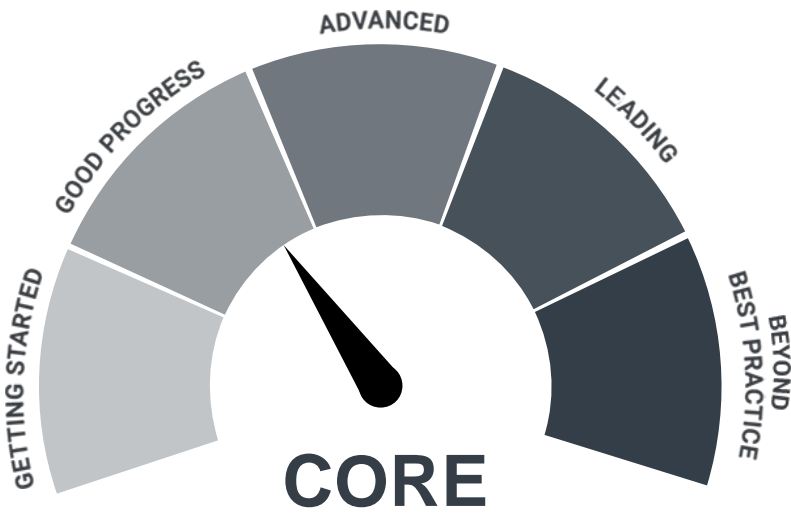


DASHBOARD



SUMMARY

For the 2020 APCO Annual Report, *Promega Corporation* has achieved Level 2 (Good Progress) for the core criteria. All seven core criteria were answered and six out of six recommended criteria were answered.

INDICATORS

This Annual Report shows the organisation's performance against the APCO Packaging Sustainability Framework.

CORE: All core criteria.

RECOMMENDED: All answered recommended criteria.

TOTAL: Sum of all core and answered recommended criteria.

LEADERSHIP: All criteria that defines leadership on packaging sustainability.

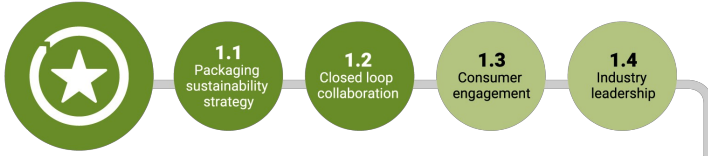
OUTCOMES: All criteria that lead to direct improvement in packaging sustainability.

OPERATIONS: All criteria related to business operations for improving packaging sustainability.

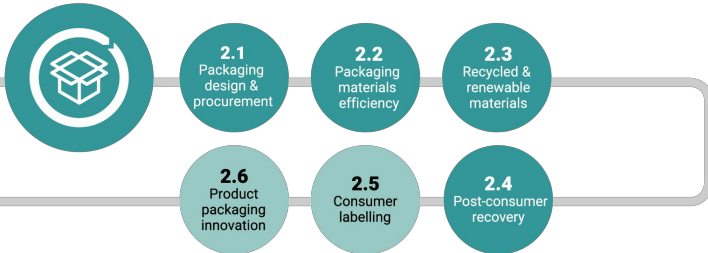
REPORTING FRAMEWORK

OVERVIEW

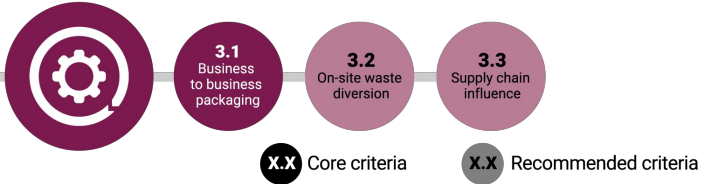
1. LEADERSHIP



2. OUTCOMES



3. OPERATIONS



LEVEL DESCRIPTION

LEVEL 0

NOT YET STARTED

The organisation does not have a plan and has not started investigating options for action under the relevant criteria.

LEVEL 1

GETTING STARTED

The organisation is developing a plan or is investigating options for action under that criteria (e.g. agreeing on goals and targets).

LEVEL 2

GOOD PROGRESS

There are documented procedures in place (e.g. to implement the Sustainable Packaging Guidelines) or some progress has been achieved, data tracking shows up to 20% of products have achieved the desired outcome.

LEVEL 3

ADVANCED

The organisation has specific, measurable targets in place, or data tracking shows that >20% of products have achieved the desired outcome.

LEVEL 4

LEADING

The organisation has more rigorous procedures in place, or an ambitious target has been met e.g. >50% of products have achieved the desired outcome.

LEVEL 5

BEYOND BEST PRACTICE

The organisation has more rigorous procedures in place, or an ambitious target has been met e.g. 100% of products have achieved the desired outcome.

ABOUT THE APCO FRAMEWORK

The APCO Packaging Sustainability Framework (the Framework) aims to provide a consistent and transparent framework for assessing and tracking packaging sustainability across organisations. The Framework consists of thirteen independent criteria that belong to three separate categories. These categories are (i) Leadership, (ii) Packaging Outcomes and (iii) Operations. The diagram above shows the different criteria included within each of these three categories.

There are two types of criteria within the Framework:

- **Core criteria** are mandatory and must be answered by all Members when completing the Annual Reporting Tool.
- **Recommended criteria** are not mandatory, but Members are encouraged to complete as many recommended criteria as possible.

Each criteria is divided into five levels of performance ranging from 'getting started' to 'beyond best practice'. The diagram above illustrates the key differences between performance levels. Please see www.packagingcovenant.org.au for more information.

ABOUT APCO

APCO is a co-regulatory, not for profit organisation that partners with government and industry to reduce the harmful impact of packaging on the Australian environment. APCO promotes sustainable design and recycling initiatives, waste to landfill reduction activities and circular economy projects. APCO Annual Reports are a key Membership obligation, providing Member organisations with a transparent method to track their packaging sustainability performance. Annual reporting also supplies APCO with valuable data and information to identify industry trends and advancements.

Promega Corporation

FREE FORM QUESTIONS

Describe initiatives, processes or practices that you have implemented during your chosen reporting period that have improved packaging sustainability.

Promega implemented changes to our kit packaging boxes that use sustainably sourced materials, reduce material use and promote recycling for customers.

CASE STUDIES

Case Study

New kit boxes featured on <https://www.promega.com/c/global/kit-packaging/> Certified as Sustainable Turn the kit box over and you will find new sustainability logos. As always, all of our paper packaging is recyclable. Now Promega uses Sustainable Forestry Initiative (SFI) Chain of Custody level certified fiber in our printed package material. The SFI label indicates a commitment to tracking fiber from forest content to end product using a third-party internationally recognized sustainability standard. In addition, the ink, coating and glue are sourced from renewable products containing no controversial materials. Our new packaging is part of our commitment to sustainability and corporate responsibility efforts that benefit our planet, people, communities and products. Less is More While our packaging materials are more sustainable, we still wanted to use less paper. By streamlining our labeling system to deliver information more efficiently, only one label is needed on the package instead of two. We re-engineered the box to be more efficiently designed, including implementing an auto-bottom configuration, meaning we can use less paperboard, ink and adhesive. Plus, the boxes will weigh less to reduce carbon emissions from shipping. Last year, Promega was recognized with the Graphic Design USA Award for Sustainable Packaging for these new kit boxes.

COMMITMENTS

The Action Plan detailed below, is a public commitment to advancing packaging sustainability outcomes as a Member of APCO. Members are encouraged to have a comprehensive internal plan with the steps they will take to fulfill their commitments.

Criteria 1.1 PACKAGING SUSTAINABILITY STRATEGY (core)

1. Getting started

Your organisation is committed to: Having a strategy to improve packaging sustainability.

Criteria 1.2 CLOSED-LOOP COLLABORATION (core)

1. Getting started

Your organisation is committed to: Investigating options to join or start a collaborative initiative to address barriers to the recovery of waste packaging.

Criteria 1.3 CONSUMER ENGAGEMENT (recommended)

4. Leading

Your organisation is committed to engaging consumers about packaging sustainability by: (1) Including packaging information on your website or other publications. (2) Applying on-pack claims or labels. (3) Designing packaging to reduce the impacts of consumption. (4) Running a marketing campaign.

Criteria 1.4 INDUSTRY LEADERSHIP (recommended)

2. Good progress

Your organisation is committed to: (1) Investigating collaborative or industry leadership actions to improve packaging sustainability. (2) Having at least one leadership initiative externally recognised through an award or other formal process.

Criteria 2.1 PACKAGING DESIGN & PROCUREMENT (core)

2. Good progress

Your organisation is committed to: Reviewing or starting to review your packaging with reference to the Sustainable Packaging Guidelines (SPG) or equivalent to identify any improvement opportunities.

COMMITMENTS

Criteria 2.2 PACKAGING MATERIALS EFFICIENCY (core)

3. Advanced

Your organisation is committed to: Reducing packaging weight or optimising packaging material efficiency in more than 20% of products.

Criteria 2.3 RECYCLED & RENEWABLE MATERIALS (core)

3. Advanced

Your organisation is committed to: Incorporating or optimising recycled and/or renewable content in more than 20% of products.

Criteria 2.4 POST-CONSUMER RECOVERY (core)

3. Advanced

Your organisation is committed to: Having more than 20% of primary packaging be recoverable through existing post-consumer recovery systems.

Criteria 2.5 CONSUMER LABELLING (recommended)

3. Advanced

Your organisation is committed to: Labelling more than 20% of products with disposal or recovery information.

Criteria 2.6 PRODUCT-PACKAGING INNOVATION (recommended)

1. Getting started

Your organisation is committed to: Investigating opportunities to change the design or delivery of at least one product to improve packaging sustainability.

Criteria 3.1 BUSINESS-TO-BUSINESS PACKAGING (core)

1. Getting started

Your organisation is committed to: Developing a plan or investigating opportunities to reduce the amount of single-use business-to-business (B2B) packaging used internally or sent to customers.

COMMITMENTS

Criteria 3.2 ON-SITE WASTE DIVERSION (recommended)

4. Leading

Your organisation is committed to: Investigating opportunities to divert more than 50% of solid waste generated at facilities (factories, warehouses, offices, retail stores etc.) from landfill.

Criteria 3.3 SUPPLY CHAIN INFLUENCE (recommended)

0. Not yet started

Your organisation has no commitment to this criteria.

SIGN OFF

Ghislaine Samways

General Manager

Wednesday, 1 July 2020

DISCLAIMER

The copyright owners of this report are The Australian Packaging Covenant Organisation (APCO) and the University of Technology Sydney (UTS). The data contained within this report is owned exclusively by APCO. This report (including all text, images and attachments) has been prepared for the exclusive use and benefit of the copyright holders and each addressee(s) only. The copyright owners give their consent and grant a limited licence to each addressee(s) for the use of this report, in its entirety, to be reproduced, distributed and communicated to third parties or made public at the discretion of the addressee(s). All other rights are strictly reserved. The addressee(s) is not permitted to reproduce, distribute or communicate separate subsections, images or individual pages of this report. The report can only be reproduced, distributed or communicated in full. The information contained in this report does not constitute investment or business advice and addressee(s) should not rely on the information for such purposes. All addressee(s) should obtain their own independent advice before making any financial decisions based on the information included in this report. 2017-2020 ©.

Suite 1102, Level 11, 55 Clarence Street, Sydney, NSW, 2000 ☎ (02) 8381 3700 ✉ apco@packagingcovenant.org.au