



75 g.m⁻²



100% Satisfaction
Product Replacement Guaranteed

5g.m⁻² Lighter
The Eco-Logical Choice

COPIER | LASER | INKJET |

ISO 9001 & ISO 14001 CERTIFIED



Herdade de Espirra Project

The **Herdade de Espirra** is a vast estate dedicated to farming and forestry activities, occupying 1,700 hectares in the municipality of Palmela, 60 km south of Lisbon, Portugal. While being located in a region that is particularly suited for vine growing, the estate is characterized by the coexistence of vineyards with other crops, fruit groves and grazing land, as well as plantations of trees such as poplars, oak trees, eucalyptuses and pine trees.

Purchased by Portucel in 1985, the Herdade de Espirra is a diversified space in terms of soil occupation. On entering the farm, the first you will see is its social area, comprising a meeting room, offices, a mansion, the cellar and the warehouse. The outline of the buildings is that of the region's traditional "monte", blending harmoniously into the landscape.

The vineyards, from which is produced the excellent Herdade de Espirra red, occupy approximately 37 hectares. The wine is made in the estate's own winery. The nurseries, which produce mainly forest species, in particular the eucalyptus, but also ornamental plants and fruit trees, have an area of 23.6 hectares. Part of the production is intended for the Group's own consumption, and the remainder is sold in the market, namely to private producers, associations and institutions.

RAIZ, the forest and paper research institute where the Portucel Soporcel group holds a majority stake, is represented at Espirra by the team responsible for the program of eucalyptus genetic improvement.

To this end, there is a nursery with modern greenhouses dedication to the cultivation of plants for testing, the development of vegetative production techniques, groves for seed production to make controlled cross-breeding, and support areas to the development of new clones.

The area of forest occupies more than 1,600 hectares, featuring a variety of farming and forestry activities, from wood and cork production, to the extraction of pine nuts, cattle grazing and apiculture. The most representative species are the eucalyptus (677 hectares), the oak tree (532 hectares), the stone pine (48 hectares), the poplar (35 hectares) and the maritime pine (2 hectares), and there are also mixed cultivation areas.

HERDADE DE ESPIRRA

Location: Palmela, Area: 1700 hectares

Soil occupation:

Vineyards (37 hectares)

Nurseries (23.6 hectares)

Eucalyptus (677 hectares)

Oak (532 hectares)

Stone pine (48 hectares)

Poplar (35 hectares)

Maritime pine (2 hectares) and mixed plantations

The Portucel Soporcel group is responsible for woodland assets up and down the country, totaling approximately 120 thousand hectares. This land is managed in accordance with the principles set out in the group's [Forestry Policy](#).

The eucalyptus, more specifically the *Eucalyptus globulus* species, which is considered at world level to provide the ideal fibre for the production of high quality papers, occupies 74% of this area.

In the [Biodiversity](#) scope, the Group has established partnerships with the WWF – World Wide Fund for Nature



The majority of the Group's forest assets are certified by the Forest Stewardship Council (FSC), whose certification system guarantees that the Company's forests are responsibly managed at environmental, economic and social level, and comply with strict and internationally recognized criteria. On a first phase the scope of FSC certification has been restricted to mainland estate, owned by or leased to the Group, and to eucalyptus round wood. Areas excluded from FSC certification comprise of some areas where long-term sustainable management can not be guaranteed by the company (some leasing contracts or areas where economic sustainability is not foreseen). Nevertheless, in these non FSC-certified areas, the Groups publicly assume it's commitment towards a responsible conduct, respecting the spirit and intention expressed in FSC's Principles and Criteria.

Besides certifying its own managed forests, the Group has been actively promoting forest certification adoption by private forest owners all over the country.

The Group's Herdade de Espirra (Espirra Estate) is a good example of its long-standing strategy focused on biodiversity and multiple soil use. It is in this estate that the Group produces the Herdade de Espirra wine, which has already won several prizes in Portugal and abroad.

Its important forest assets make the Group a net carbon fixer. By the end of 2007, the plantations managed by the Company had an accumulated retention of 7.7 million tons of CO₂.

The Group sees the forest as a strategic source of wealth for the country: in fact the national forest based industry generates 2.8% of the Portuguese Gross Domestic Product (GDP).

The Portucel Soporcel group is strongly committed to applied research, an activity developed at [RAIZ](#) , Instituto de Investigação da Floresta e Papel (Forest and Paper Research Institute). [RAIZ](#)  supports the forestry and paper industries, providing technological support and specialized training.

Eco-Logical | 75 g.m⁻²

MULTIFUNCTIONAL

5g.m⁻² Lighter, The Eco-Logical Choice.

THE ECOLOGICAL CHOICE

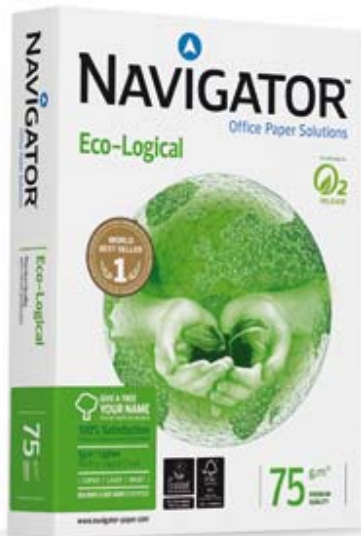
Uses fewer natural resources whilst guaranteeing a premium quality performance.

GIVE A TREE YOUR NAME

The opportunity to give your name to one of the several million trees grown every year in our forest nurseries.

100% SATISFACTION

High quality, cost efficient, office paper backed by a Product Replacement Guarantee.



PRESERVING BIO-DIVERSITY



CONTRIBUTING TOWARDS PRESERVING THE ENVIRONMENT

grupo Portucel Soporcel integrates biodiversity conservation in its forest management model, built upon partnerships with environmental organizations.



By buying products with an FSC label you are supporting the growth of responsible forest management worldwide.

GREEN INNOVATION



FIGHTING CLIMATE CHANGE:

Our forests store the equivalent to the CO₂ emissions generated by 1,5 million cars driving around the Globe, and each year fixed more than twice the fossil CO₂ produced in all our mills.

Our Give a Tree Your Name program not only promotes responsible reforestation, giving you a real part in the prevention of global warming, but also contributes to biodiversity protection and preservation.

Increased usage of renewable energy. More than 75% of the total energy produced by grupo Portucel Soporcel is obtained from renewable resources.

Eco-Logical

75g.m⁻²

A4

210x297 mm
500 SHEETS / REAM | 5 REAMS / BOX

A3

420x297 mm
500 SHEETS / REAM | 5 REAMS / BOX

A4

FastPack
2500 SHEETS
Unwrapped



75 g.m⁻²



**GIVE A TREE
YOUR NAME**

100% Satisfaction
Product Replacement Guaranteed

5g.m⁻² Lighter
The Eco-Logical Choice

| **COPIER** | **LASER** | **INKJET** |

ISO 9001 & ISO 14001 CERTIFIED

[HOME](#)
[KEY FEATURES](#)
[RANGE](#)
[WHERE TO BUY](#)
[MEDIA CENTRE](#)
[CONTACT US](#)